

TMT10plex Mass Tag Labeling Kits and Reagents

Pub. No. MAN0016969

Rev B.0

Pub. Part No. 2162457.5

Number	Description
90110	TMT10plex Isobaric Label Reagent Set , sufficient reagents for one 10plex isobaric experiment Contents: TMT¹⁰-126TM Label Reagent , 1 × 0.8mg TMT¹⁰-127NTM Label Reagent , 1 × 0.8mg TMT¹⁰-127CTM Label Reagent , 1 × 0.8mg TMT¹⁰-128NTM Label Reagent , 1 × 0.8mg TMT¹⁰-128CTM Label Reagent , 1 × 0.8mg TMT¹⁰-129NTM Label Reagent , 1 × 0.8mg TMT¹⁰-129CTM Label Reagent , 1 × 0.8mg TMT¹⁰-130NTM Label Reagent , 1 × 0.8mg TMT¹⁰-130CTM Label Reagent , 1 × 0.8mg TMT¹⁰-131TM Label Reagent , 1 × 0.8mg
90111	TMT10plex Isobaric Label Reagent Set , sufficient reagents for three 10plex isobaric experiments Contents: TMT¹⁰-126 Label Reagent , 3 × 0.8mg TMT¹⁰-127N Label Reagent , 3 × 0.8mg TMT¹⁰-127C Label Reagent , 3 × 0.8mg TMT¹⁰-128N Label Reagent , 3 × 0.8mg TMT¹⁰-128C Label Reagent , 3 × 0.8mg TMT¹⁰-129N Label Reagent , 3 × 0.8mg TMT¹⁰-129C Label Reagent , 3 × 0.8mg TMT¹⁰-130N Label Reagent , 3 × 0.8mg TMT¹⁰-130C Label Reagent , 3 × 0.8mg TMT¹⁰-131 Label Reagent , 3 × 0.8mg
90406	TMT10plex Isobaric Label Reagent Set , sufficient reagents for one 10plex isobaric experiment Contents: TMT¹⁰-126 Label Reagent , 1 × 5mg TMT¹⁰-127N Label Reagent , 1 × 5mg TMT¹⁰-127C Label Reagent , 1 × 5mg TMT¹⁰-128N Label Reagent , 1 × 5mg TMT¹⁰-128C Label Reagent , 1 × 5mg TMT¹⁰-129N Label Reagent , 1 × 5mg TMT¹⁰-129C Label Reagent , 1 × 5mg TMT¹⁰-130N Label Reagent , 1 × 5mg TMT¹⁰-130C Label Reagent , 1 × 5mg TMT¹⁰-131 Label Reagent , 1 × 5mg

-
- 90113** **TMT10plex Isobaric Mass Tag Labeling Kit**, sufficient reagents for three 10plex isobaric experiments
- Contents:**
- TMT¹⁰-126 Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-127N Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-127C Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-128N Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-128C Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-129N Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-129C Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-130N Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-130C Label Reagent**, 3 × 0.8mg
 - TMT¹⁰-131 Label Reagent**, 3 × 0.8mg
 - Dissolution Buffer** (1M triethyl ammonium bicarbonate), 5mL
 - Denaturing Reagent** (10% SDS), 1mL
 - Reducing Reagent** (0.5M TCEP), 1mL
 - Iodoacetamide**, 12 × 9mg
 - Quenching Reagent** (50% hydroxylamine), 1mL
 - Pierce Trypsin Protease, MS Grade**, 5 × 20µg
 - Trypsin Storage Solution**, 250µL
 - Albumin, Bovine**, 2.5mg
- A34807** **TMT11-131C**, 1 × 5mg
- A34808** **TMT10plex Isobaric Label Reagent Set plus TMT11-131C**, sufficient reagents for one 11plex isobaric experiment
- Contents:**
- TMT¹⁰-126 Label Reagent**, 1 × 5mg
 - TMT¹⁰-127N Label Reagent**, 1 × 5mg
 - TMT¹⁰-127C Label Reagent**, 1 × 5mg
 - TMT¹⁰-128N Label Reagent**, 1 × 5mg
 - TMT¹⁰-128C Label Reagent**, 1 × 5mg
 - TMT¹⁰-129N Label Reagent**, 1 × 5mg
 - TMT¹⁰-129C Label Reagent**, 1 × 5mg
 - TMT¹⁰-130N Label Reagent**, 1 × 5mg
 - TMT¹⁰-130C Label Reagent**, 1 × 5mg
 - TMT¹⁰-131 Label Reagent**, 1 × 5mg
 - TMT11-131C Label Reagent**, 1 × 5mg
- A37724** **TMT11-131C**, 3 × 0.8mg

A37725 TMT10plex Isobaric Label Reagent Set plus TMT11-131C, sufficient reagents for three 11plex isobaric experiments

Contents:

- TMT¹⁰-126 Label Reagent, 3 × 0.8mg
- TMT¹⁰-127N Label Reagent, 3 × 0.8mg
- TMT¹⁰-127C Label Reagent, 3 × 0.8mg
- TMT¹⁰-128N Label Reagent, 3 × 0.8mg
- TMT¹⁰-128C Label Reagent, 3 × 0.8mg
- TMT¹⁰-129N Label Reagent, 3 × 0.8mg
- TMT¹⁰-129C Label Reagent, 3 × 0.8mg
- TMT¹⁰-130N Label Reagent, 3 × 0.8mg
- TMT¹⁰-130C Label Reagent, 3 × 0.8mg
- TMT¹⁰-131 Label Reagent, 3 × 0.8mg
- TMT11-131C Label Reagent, 3 × 0.8mg

Storage: Upon receipt store at -20°C. Reagents are shipped with dry ice.

Note: Products are for research use only – do not use for diagnostic procedures.

Table of Contents

Table of Contents.....	3
Introduction	4
Procedure Summary.....	4
Important Product Information	5
Additional Materials Required.....	6
Material Preparation	6
Preparing and Labeling Peptides with the TMT Isobaric Mass Tags	6
Troubleshooting.....	8
Additional Information	8
A. Data Acquisition Methods	8
B. Data Analysis and Quantitation	9
Related Thermo Scientific Products	9
General References	10
Limited Use Label License: TMT10plex™ Isobaric Mass Tag Labeling Kits and Reagents.....	10

Introduction

The Thermo Scientific™ TMT™ Mass Tag Labeling Kits and Reagents enable multiplex relative quantitation by mass spectrometry (MS). Each mass-tagging reagent within a set has the same nominal mass (i.e., isobaric) and chemical structure composed of an amine-reactive NHS-ester group, a spacer arm and a mass reporter (Figure 2A). The reagent set can be used to label up to 11 different peptide samples prepared from cells or tissues. For each sample, a unique reporter mass (i.e., 126-131Da) in the low mass region of the MS/MS spectrum is used to measure relative protein expression levels during peptide fragmentation (Figure 1).

The Thermo Scientific™ TMT10plex™ Label Reagents and TMT11-131C share an identical structure with Thermo Scientific™ TMTzero™ and TMTsixplex™ Reagents, but contain different numbers and combinations of ¹³C and ¹⁵N isotopes in the mass reporter. The different isotopes result in a set of tags that have 6 mDa monoisotopic mass differences in the reporter that can be detected using high resolution Thermo Scientific™ Orbitrap™ Mass Spectrometry Instruments. Advantages of the TMT Label Reagents include increased sample multiplexing for relative quantitation, increased sample throughput and fewer missing quantitative channels among samples.

Procedure Summary

Protein extracts isolated from cells or tissues are reduced, alkylated and digested overnight. Samples are labeled with the TMT Reagents and then mixed before sample fractionation and clean-up. Labeled samples are analyzed by high resolution Orbitrap LC-MS/MS before data analysis to identify peptides and quantify reporter ion relative abundance (Figure 2).

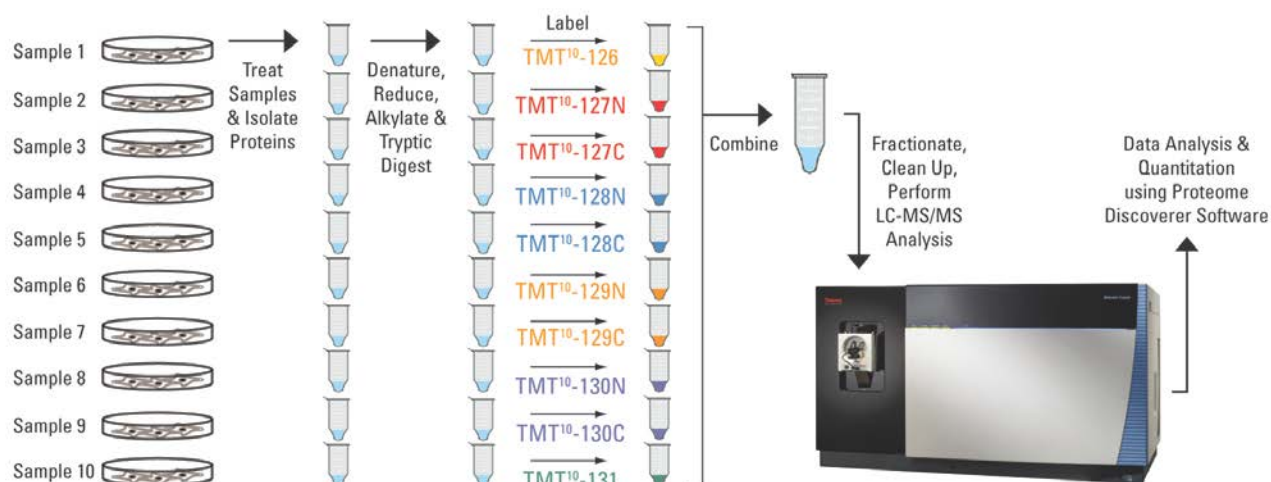


Figure 1. Procedure schematic for using the Thermo Scientific TMT10plex Label Reagents.

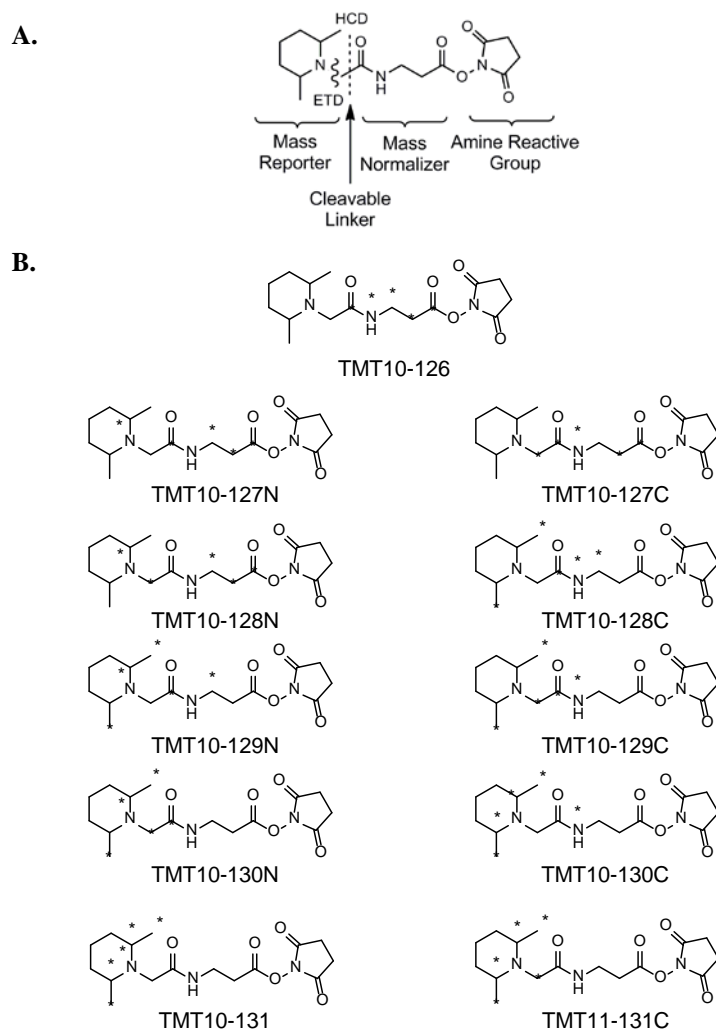


Figure 2. Chemical structures of the Thermo Scientific TMT Label Reagents.

A. Functional regions of the reagent structure including MS/MS fragmentation sites by higher-energy collision dissociation (HCD) and electron transfer dissociation (ETD).

B. TMT10plex reagent and TMT11-131C reagent structures and isotope positions (*).

Important Product Information

- The TMT reagents are highly moisture-sensitive. To avoid moisture condensation onto the product, the vial must be equilibrated to room temperature before opening. Store unused reagent in foil pouch with desiccant at -20°C.
- The TMT reagents are amine-reactive and modify lysine residues and peptide N-termini. All amine-containing buffers and additives must be removed before digestion and labeling.
- All samples must be digested, labeled and then mixed equally before desalting, fractionation and LC-MS/MS. For optimal results, use 25-100µg of peptide for each labeling reaction.
- To avoid contamination of MS samples, always wear gloves when handling samples and gels. Use ultrapure MS-grade reagents. Perform sample preparation in a clean work area.
- The TMTzero Label Reagent (Product No. 90067) can be used to optimize methods before multiplexed analysis of samples with TMT10plex Label Reagent sets.

Additional Materials Required

- Microcentrifuge tubes
- Anhydrous acetonitrile (Acetonitrile, LC-MS Grade, Product No. 51101)
- Water, LC-MS Grade (Product No. 51140)
- Chilled (-20°C) acetone
- Protein assay (e.g., Thermo Scientific™ BCA Protein Assay Kit, Product No. 22235)
- 75-300µm capillary C₁₈ reversed-phase column
- High-resolution Orbitrap Mass Spectrometer with online liquid chromatography system (see Table 1 for recommended instruments)
- Data analysis software (e.g., Thermo Scientific™ Proteome Discoverer™ Software)
- Optional: C18 spin tips or columns (e.g., Thermo Scientific™ Pierce™ C18 Spin Columns, Product No. 89870 or Pierce™ C18 Tips, Product No. 87784)

Material Preparation

Note: The 50% hydroxylamine and 10% SDS stock solutions provided with the kit may precipitate during storage. Warm both solutions to room temperature and vortex before use. The amounts listed below are sufficient for preparing and labeling 10 samples.

100mM TEAB (triethyl ammonium bicarbonate)	Add 500µL of the Dissolution Buffer (1M TEAB) to 4.5mL of ultrapure water.
Lysis Buffer	Add 200µL of the Denaturing Reagent (10% SDS) to 1.8mL of 100mM TEAB.
200mM TCEP	Add 70µL of the Reducing Reagent (0.5M TCEP) to 70µL of ultrapure water. Then add 35µL of the Dissolution Buffer (1M TEAB).
5% Hydroxylamine	Add 50µL of the Quenching Reagent (50% hydroxylamine) to 450µL of 100mM TEAB.

Preparing and Labeling Peptides with the TMT Isobaric Mass Tags

Note: BSA can be used as a control sample for method optimization. Dissolve BSA to 1mg/mL using 100mM TEAB, pH 8.5. HEPES pH 8.5 buffer can be used as alternative buffer for digestion and labeling; however, C18 desalting is required to remove the buffer. Use 25-100µg of protein per labeling reaction. The Thermo Scientific™ Pierce™ Mass Spec Sample Prep Kit for Cultured Cells can also be used to prepare peptide digests for TMT reagent labeling.

A. Preparing Whole Cell Protein Extracts

1. Culture cells to harvest at least 100µg of protein per condition. For best results, culture a minimum of 2×10^6 cells.
Note: Rinse cells 2-3 times with 1X PBS to remove cell culture media. Pellet cells using low-speed centrifugation (i.e., $< 1000 \times g$) to prevent premature cell lysis.
2. Lyse the cells by adding five cell-pellet volumes of Lysis Buffer (i.e., 100µL of Lysis Buffer for a 20µL cell pellet).
Note: Lysis buffers such as 8M urea (Product No. 29700) in 50mM TEAB or HEPES buffer, pH 8 may be used as alternative denaturing cell lysis buffers. For urea-based lysis buffer, protein samples must be diluted to $< 1M$ urea before digestion, and the final C18 desalting step (C.6) is not optional. Addition of protease and/or phosphatase inhibitors during lysis is optional and may interfere with MS analysis.
Note: Depending on the Lysis Buffer used it may be necessary to reduce sample viscosity by shearing DNA using a microtip sonicator or addition of a nuclease (e.g., Thermo Scientific™ Pierce™ Universal Nuclease for Cell Lysis, Product No. 88700)
3. Centrifuge lysate at $16,000 \times g$ for 10 minutes at 4°C.
4. Carefully separate the supernatant and transfer into a new tube.

-
- Determine the protein concentration of the supernatant using established methods such as the BCA Protein Assay Kit (Product No. 23227).

Note: Use samples at ≥ 2 mg/mL. Less concentrated samples may be used; however, it might be necessary to use larger volumes of reducing/alkylating reagents.

- Transfer 100 μ g per condition into a new tube and adjust to a final volume of 100 μ L with 100mM TEAB.
- Add 5 μ L of the 200mM TCEP and incubate sample at 55°C for 1 hour.
- Immediately before use, dissolve one tube of iodoacetamide (9mg) with 132 μ L of 100mM TEAB to make 375mM iodoacetamide. Protect solution from light.
- Add 5 μ L of the 375mM iodoacetamide to the sample and incubate for 30 minutes protected from light at room temperature.
- Add six volumes (~ 600 μ L) of pre-chilled (-20°C) acetone. Allow the precipitation to proceed overnight.
Note: Methanol/chloroform is the recommended solvent for precipitation of proteins derived from tissue extracts.
- Centrifuge the samples at 8000 $\times g$ for 10 minutes at 4°C. Carefully invert the tubes to decant the acetone without disturbing the white pellet. Allow the pellet to dry for 2-3 minutes.

B. Protein Digestion

- Resuspend 100 μ g of acetone-precipitated (or lyophilized) protein pellets with 100 μ L of 100mM TEAB or HEPES pH 8.5.
Note: An acetone-precipitated pellet may not completely dissolve; however, after proteolysis at 37°C, all the protein (peptides) will be solubilized.
- Immediately before use, add 20 μ L of the Trypsin Storage Solution to the bottom of the trypsin glass vial and incubate for 5 minutes. Store any remaining reagent in single-use volumes at -80°C (e.g., 2.5 μ g of trypsin per 100 μ g of protein).
- Add 2.5 μ L of trypsin (i.e., 2.5 μ g) per 100 μ g of protein. Digest the sample overnight at 37°C.

C. Peptide Labeling

- Immediately before use, equilibrate the TMT Label Reagents to room temperature. For the 0.8mg vials, add 41 μ L of anhydrous acetonitrile to each tube. For the 5mg vials, add 256 μ L of solvent to each tube. Allow the reagent to dissolve for 5 minutes with occasional vortexing. Briefly centrifuge the tube to gather the solution.
Note: Reagents dissolved in anhydrous acetonitrile are stable for one week when stored at -20°C and warmed to room temperature before opening. Anhydrous ethanol can be used as an alternative solvent to dissolve reagents.
- Optional: Measure protein digest concentration using Thermo Scientific™ Pierce™ Quantitative Fluorescent Peptide Assay (Product No. 23290) or Thermo Scientific™ Pierce™ Quantitative Colorimetric Peptide Assay (Product No. 23275).
- Carefully add 41 μ L of the TMT Label Reagent to each 100 μ L sample (25-100 μ g protein digest). Alternatively, transfer the reduced and alkylated protein digest to the TMT Reagent vial.
Note: Labeling more than 100 μ g of protein digest per reaction requires additional TMT Label Reagent.
- Incubate the reaction for 1 hour at room temperature.
- Add 8 μ L of 5% hydroxylamine to the sample and incubate for 15 minutes to quench the reaction.
- Combine equal amounts of each sample in a new microcentrifuge tube and speedvac to dry labeled peptide sample.
- Clean-up samples using peptide desalting columns (Product No. 89852) or equivalent before high-resolution LC-MS analysis. Alternatively, Thermo Scientific™ Pierce™ High pH Reversed-Phase Peptide Fractionation Kit (Product No. 84868) can be used to clean up and fractionate TMT-labeled peptides to increase the number of peptide identifications.
Note: TMT-labeled peptide concentration can be measured after clean up using the Pierce Quantitative Colorimetric Peptide Assay. The Pierce Quantitative Fluorescent Peptide Assay cannot be used to measure TMT-labeled peptide concentrations.

Troubleshooting

Problem	Possible Cause	Solution
Poor labeling	A primary amine-based buffer was used (e.g., Tris, glycine)	Use non-primary amine-based buffers (e.g., TEAB, HEPES)
	Incorrect buffer pH	Make sure the buffer pH is ~8.0-8.5
	Too much sample was used	Label 25-100µg sample per 0.8mg of TMT reagent
	Incorrect solvent was used	Use dry acetonitrile or ethanol to reconstitute tags
	Reagents hydrolyzed	Avoid exposing tags to moisture Store unused reagents with desiccant at -20°C
Poor protein quantitation	Incorrect instrument method used	Optimize TMT reporter ion MS/MS fragmentation
	Too little sample analyzed	Increase sample amount and optimize ion injection
	Poor chromatography	Optimize LC gradient to maximize MS/MS of unique peptides
	Co-isolation of peptides during MS	Reduce sample complexity by pre-fractionating peptides
		Decrease quadrupole isolation width if applicable
	Use MS3 methods (i.e. SPS-MS3)	

Additional Information

A. Data Acquisition Methods

Quantitation of peptides labeled with Thermo Scientific™ Tandem Mass Tag™ Reagents requires a high-resolution Orbitrap Mass Spectrometer capable of MS/MS fragmentation (Table 1). To resolve near-isobaric reporter ions, MS/MS resolution must be > 50,000 at 150 *m/z*. Higher-energy collision dissociation (HCD) is recommended for TMT10plex reporter ion fragmentation. Optimal HCD fragmentation energy is instrument-dependent and can be optimized using TMTzero Reagents. Electron transfer dissociation (ETD) may be used as an alternative fragmentation method for peptide identification and quantitation; however, ETD is not recommended for TMT10plex Reagents because of reporter ion overlap (Table 2).

Table 1. Instruments and MS/MS fragmentation options for peptide identification and quantitation with Thermo Scientific TMT Reagents.

<u>Instrument</u>	<u>Fragmentation Method</u>	<u>Minimum Resolution Setting</u>	<u>Reference(s)</u>
Thermo Scientific Orbitrap Fusion™ Tribrid™ Mass Spectrometer	HCD/SPS-MS3	60,000	McAllister, <i>et al.</i> (2014)
Thermo Scientific Orbitrap Elite™ Mass Spectrometer	HCD/MS3	30,000	Viner, <i>et al.</i> (2012)
Thermo Scientific Q Exactive™ Mass Spectrometer	HCD/MS2	35,000	Wühr, <i>et al.</i> (2012)
Thermo Scientific Orbitrap Velos Pro Mass Spectrometer	HCD/MS2	30,000	Ting, <i>et al.</i> (2011), Wenger, <i>et al.</i> (2011)

B. Data Analysis and Quantitation

The peptide mass modification by the TMT10plex Reagents is identical to TMTsixplex Reagents and present in the UNIMOD database (www.unimod.org) and are listed below. Proteome Discoverer Software (2.1 and above) is recommended for TMT10plex relative quantitation. Additional software programs that may be used for TMT quantitation include Matrix Science™ Mascot™ Software (2.5 and above) and Proteome Software™ Scaffold™ Q+ Software. For data acquired using a combination of fragmentation methods (i.e., HCD/MS3 or HCD/ETD), Proteome Discoverer Software may be necessary to merge search results.

Table 2. Modification masses of the Thermo Scientific TMT Label Reagents.

<u>Label Reagent</u>	<u>Label Reagent</u>	<u>Modification Mass (monoisotopic)</u>	<u>Modification Mass (average)</u>	<u>HCD Monoisotopic Reporter Mass*</u>	<u>ETD Monoisotopic Reporter Mass**</u>
TMT ¹⁰ -126	TMT ⁶ -126	229.162932	229.2634	126.127726	114.127725
TMT ¹⁰ -127N	TMT ⁶ -127	229.162932	229.2634	127.124761	115.124760
TMT ¹⁰ -127C	-	229.162932	229.2634	127.131081	114.127725
TMT ¹⁰ -128N	-	229.162932	229.2634	128.128116	115.124760
TMT ¹⁰ -128C	TMT ⁶ -128	229.162932	229.2634	128.134436	116.134433
TMT ¹⁰ -129N	TMT ⁶ -129	229.162932	229.2634	129.131471	117.131468
TMT ¹⁰ -129C	-	229.162932	229.2634	129.137790	116.134433
TMT ¹⁰ -130N	-	229.162932	229.2634	130.134825	117.131468
TMT ¹⁰ -130C	TMT ⁶ -130	229.162932	229.2634	130.141145	118.141141
TMT ¹⁰ -131	TMT ⁶ -131	229.162932	229.2634	131.138180	119.138176
TMT11-131C		229.169252	229.2634	131.144499	118.141141

* HCD is a collisional fragmentation method that generates ten unique reporter ions from 126 to 131Da.

**ETD is a non-ergodic fragmentation method that generates six unique reporter ions from 114 to 119Da.

Related Thermo Scientific Products

90114	1M Triethylammonium bicarbonate (TEAB), 50mL
90115	50% Hydroxylamine, 5mL
90067	TMTzero Label Reagent, 5 × 0.8mg
90061	TMTsixplex Isobaric Label Reagent Set, 1 × 0.8mg
90064	TMTsixplex Isobaric Mass Tagging Kit
90308	TMTsixplex Isobaric Label Reagent Set, 16 × 0.2mg
90309	TMT10plex Isobaric Label Reagent Set, 8 × 0.2mg
90100	iodoTMTzero™ Label Reagent, 5 × 0.2mg
90101	iodoTMTsixplex™ Label Reagent Set, 1 × 0.2mg
90103	iodoTMTsixplex Isobaric Mass Tag Labeling Kit
84840	Pierce™ Mass Spec Sample Prep Kit for Cultured Cells
23275	Pierce Quantitative Fluorescent Peptide Assay
23290	Pierce Quantitative Colorimetric Peptide Assay
90057	Pierce Trypsin Protease, MS Grade
90051	Lys-C Protease, MS Grade
A32992	High Select™ Fe-NTA Phosphopeptide Enrichment Kit
A32993	High Select™ TiO ₂ Phosphopeptide Enrichment Kit
84868	Pierce High pH Reversed-Phase Peptide Fractionation Kit
89852	Pierce Peptide Desalting Columns, 25 columns
28904	Trifluoroacetic Acid, Sequanal Grade

General References

- Altelaar A.F., *et al.* (2013). Benchmarking stable isotope labeling based quantitative proteomics. *J Proteomics* **88**:14-26.
- Dillon, R., *et al.* (2011). Discovery of a novel B-Raf fusion protein related to c-Met drug resistance. *J Proteome Res* **10**(11):5084-94.
- Erikson, B.K., *et al.* (2015). Evaluating multiplexed quantitative phosphopeptide analysis on a hybrid quadrupole mass filter/linear ion trap/orbitrap mass spectrometer. *Anal Chem* **87**(2):1241-9.
- Keshishian, H., *et al.* (2015). Multiplexed, quantitative workflow for sensitive biomarker discovery in plasma yields novel candidates for early myocardial injury. *Mol Cell Proteomics*. 2015 Feb 27. pii: mcp.M114.046813
- Lee, M.V., *et al.* (2011). A dynamic model of proteome changes reveals new roles for transcript alteration in yeast. *Mol Syst Biol* **7**: 514.
- McAllister, G.C., *et al.* (2014). MultiNotch MS3 enables accurate, sensitive, and multiplexed detection of differential expression across cancer cell line proteomes. *Anal Chem* **86**(14):7150-8.
- McAllister, G.C., *et al.* (2012). Increasing the multiplexing capacity of TMTs using reporter ion isotopologues with isobaric masses. *Anal Chem* **84**(17):7469-78.
- Murphy, J.P., *et al.* (2014). Combining amine metabolomics and quantitative proteomics of cancer cells using derivatization with isobaric tags. *Proteomics* **86**(7):3585-93.
- Paulo, J.A., *et al.* (2014). A comprehensive proteomic and phosphoproteomic analysis of yeast deletion mutants of 14-3-3 orthologs and associated effects of rapamycin. *Nature* (2-3):474-86.
- Savitski, M.M., *et al.* (2014). Tracking cancer drugs in living cells by thermal profiling of the proteome. *Science* **346**(6205):1255784
- Ting, L., *et al.* (2011). MS3 eliminates ratio distortion in isobaric multiplexed quantitative proteomics. *Nat Methods* **8**: 937-40.
- Wenger, C.D., *et al.* (2011). Gas-phase purification enables accurate, multiplexed proteome quantification with isobaric tagging. *Nat Methods* **8**(11):933-5.
- Werner, T., *et al.* (2012). High-resolution enabled TMT8-plexing. *Anal Chem* **84**(16):7188-94.
- Wu, L., *et al.* (2013). Variation and genetic control of protein abundance in humans. *Nature* **499**(7456):79-82
- Wühr, M., *et al.* (2012). Accurate multiplexed proteomics at the MS2 level using the complement reporter ion cluster. *Anal Chem* **84**(21):9214-21.
- Viner, R.I., *et al.* (2013). Increasing the multiplexing of protein quantitation from 6- to 10-Plex with reporter ion isotopologues. PN_ASMS_W617_RViner_R1.
- Viner, R.I., *et al.* (2012). Relative quantitation of TMT-labeled proteomes – Focus on sensitivity and precision. Application note #566.
- Viner, R.I., *et al.* (2009). Quantification of post-translationally modified peptides of bovine α -crystallin using tandem mass tags and electron transfer dissociation. *J Proteomics* **72**(5):874-85.
- Xiong, L., *et al.* (2011). Mass spectrometric studies on epigenetic interaction networks in cell differentiation. *J Biol Chem* **286**(15):13657-68.
- Zhang, T., *et al.* (2010). Improving quantitation of TMT-labeled peptides using stepped higher-energy collisional dissociation. Application note #483.

Limited Use Label License: TMT10plex™ Isobaric Mass Tag Labeling Kits and Reagents

(Thermo Scientific Product Nos. 90110, 90111, 90113, 90406, A34807, and A34808).

The purchase of this product conveys to the buyer the non-exclusive, non-transferable right to use the product and components of the product for internal research purposes of the buyer only, whether the buyer is an academic or for-profit entity. This product is not available for use in diagnostic procedures.

The buyer must obtain a license for use of the product in performing commercial services or for any other commercial purposes. For information on purchasing a license, contact proteomics.licensing@thermofisher.com, Pierce Biotechnology, Inc., 3747 North Meridian Road, Rockford, IL 61101.

If the buyer is not willing to accept the limitations of this limited use statement, Pierce Biotechnology, Inc. is willing to accept return of the unused product with a full refund. By opening and using this product, the buyer agrees to be bound by the conditions of this Limited Use Label License.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. Tandem Mass Tag, TMT, TMTzero, TMTsixplex, TMT10plex, iodoTMTzero and iodoTMTsixplex are trademarks of Proteome Sciences plc. Unless otherwise indicated, all other trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.