Quantitative Analysis of Large Phosphopeptide Datasets Using Proteome Discoverer 2.0

David M. Horn,¹ Ilyas Singec,² Laurence Brill² ¹Thermo Fisher Scientific, San Jose, CA, USA ; ²Sanford-Burnham Research Institute, La Jolla, CA

а

Overview

Purpose: Interpretation of a complex quantitative phosphoproteomic dataset in the Proteome Discoverer 2.0 platform.

Methods: Phosphopeptides were enriched using IMAC or TiO2 from hESC and hNSC samples and the flow through, wash and enriched fractions were collected. All MS data were acquired on a Thermo Scientific[™] LTQ Orbitrap[™] Mass Spectrometer system equipped with ETD. Data were analyzed using the new study management features in Thermo Scientific[™] Proteome Discoverer[™] software 2.0 using a label-free quantification approach

Results: Proteome Discoverer 2.0 software processed the 4 stem cell samples in ~7 days with identification and quantification of phosphoproteins and phosphopeptides. Several proteins and phosphopeptides are shown to be differentially expressed and these are known to be regulated in stem cells as well as some novel proteins not known to be differentially regulated.

Introduction

With advances in liquid chromatography/mass spectrometry (LC/MS) technology, The complexity of proteomics data is increasing rapidly. It is becoming increasingly common to find datasets with 100's of GB of raw data files for complex scientific studies, putting an increased burden on downstream software tools for interpretation of such datasets.

The latest release in the Proteome Discoverer platform has several new features for analysis of complex datasets. The first major feature is the new Consensus workflow that creates persistent reports that open very large datasets quickly. Secondly, the results are presented in a new hierarchical format with linked views for protein groups, proteins, peptides, and peptide spectrum matches (PSMs). Third, the most critical feature for analysis of large quantitative datasets is the new study management. These will be demonstrated on a large dataset of a quantitative comparison between human embryonic stem cell (hESC) and neural stem cell (hNSC) derivatives

Methods

Proteins were extracted from hESCs and hNSCs, reduced and alkylated using iodoacetamide, digested using trypsin, and separated into 32 fractions using strong cation exchange (SCX) chromatography. Phosphopeptides enrichment was preformed, also collecting the flow through and wash fractions from the SCX fractions. More details on the cells and the sample preparation will be available in a forthcoming publication. Each of the fractions were analyzed in duplicate using a data dependent decision tree LC/MS/MS method on an LTQ Orbitrap Velos mass spectrometer equipped with ETD.

These study factors will be used by Proteome Discoverer to determine which quantification values will be calculated and shown in the final report.

				Sample Identifier		inple Type				Enrichme		Enrichment ste		
dy Factors			8	•	- 6		•		•	•	_	Θ ·	Θ	
				3665_TK02_0_1		ample		3955		T/02		Enriched ·		
Enrichment	Edit 🗙	8.5		3955_TIO2_e_2		angle		3955		TO2		Enriched ·		
Enrichment	Eolt X	85		3955_TK02_bv_1		ample	•	3955		TO2		Flow throug .		
	IMAC	*5		3955_TIO2_tw_2		emple		3955		102		Fiber Broug +		
	TiO2	8.5		3955_MAC_A_1		ample		3955		MAC		Enriched ·		
)	85	3	3955_M4C_e_2	ş	emple		3965		INVC		Enriched ·	2	
		85		3955_M4C_hr_1	S	ample		3055		M4C		Flow throug +	1	
nrichment step	Edit x	85	8	3955_MAC_bv_2	s	ingle		3955		MAC		Flow throug +	2	
		85	8	3956_M4C_e_1	5	smple		3956		IN4C		Enriched ·	1	
	Enriched	85	06	3956_M4C_e_2	S	engle		3956		MAC		Enriched ·	z	
F	low through/wash	85	11	3956_M4C_thr_1	S	mple		3956		M4C		Fine throug +	1	
		85	12	3956_M4C_br_2	s	enple		3956		IN/C		Flow throug +	2	
	Edit ×	85	13	3956_TK02_bv_1	s	ample		3956		T/02		Flow throug -	1	
mple name		85	14	3956_TiO2_thv_2	s	angle		3956		T/02		Flow Broug .	2	-
			15	3956_TIC2_e_1	5	engle		3956		TO2		Enriched ·	1	
	3956	8.5	76	3996_TIQ2_e_2	S	ample		3956		102		Ennched ·	2	
	3957		17	3557_102_e_1	s	mple		3957		TO2		Enriched •	1	
	3958	* 5	18	3957_TIO2_+_2	s	engle		3957		TO2		Enriched -	2	
		85	19	3957_TK02_thv_1	s	mple		3957		T/02		Flow throut .	1	
			20	3957_TIQ2_Bw.2	5	angle		3957		102		Flow throug +	2	
chnical Replicate	Edit 🗙		21	3958_TIC2_e_1	s	mple		3958		TO2		Enriched ·		
		*5		3958 TIO2 e 2	s	angle		3958		102		Ernshed ·	2	
	1			3558 TK02 fby 1	S	mole		3958		102		Fine these +	1	
	2	* 5		3958_TIO2_bv_2		male		3958	-	TO2		Flow throus -		-

FIGURE 1a) Study factors created for this analysis. b) List of samples imported into Proteome Discoverer software and the study factors assigned to those samples. Each of these samples corresponds to at least 25 raw data files.

The next step is to choose the raw files. Proteome Discoverer 2.0 software includes a new "Add Fractions" feature that groups a set of raw data files as a single sample. For this study, each set of raw data files that fit the criteria of the study factors above such as Sample 3955, TiO2-enriched phosphopeptides, technical replicate 1 are loaded simultaneously as a single sample. A typical group has 25-50 data files, with the larger collection of data files corresponding to samples where the flow through and wash samples from phoshpopeptide enrichment were run as separate LC/MS/MS runs. Once the datasets were imported into the software, the study factors defined in Figure 1 a were assigned to each sample as shown in Figure 1b above.



The next step is to choose the raw files. Proteome Discoverer 2.0 software includes a new "Add Fractions" feature that groups a set of raw data files as a single sample. For this study, each set of raw data files that fit the criteria of the study factors above such as Sample 3955, TiO2-enriched phosphopeptides, technical replicate 1 are loaded simultaneously as a single sample. A typical group has 25-50 data files, with the larger collection of data files corresponding to samples where the flow through and wash samples from phoshpopeptide enrichment were run as separate LC/MS/MS runs. Once the datasets were imported into the software, the study factors defined in Figure 1a were assigned to each sample as shown in Figure 1b above.

The third step is to create a new analysis, which includes two node-based workflows as well as the selection of the quantification details. Proteome Discoverer 2.0 software introduces a new dual workflow setup that includes a Processing Workflow for peptide identification, peptide quantification, FDR calculation and PTM site localization and adds a new Consensus workflow that is used to perform protein inference, filter the results based on FDR and other calculations, and summarize quantification results. The two workflows used for this study are shown in Figures 2a-b below.



FIGURE 2. Processing and consensus workflows for the phosphopeptide analysis.

The processing workflow in Figure 2 includes Sequest HT nodes to interpret both CID and ETD MS/MS and includes the Event Detector and Precursor Ions Area Detector nodes to calculate peak areas for identified peptides. All Sequest HT searches used 10 ppm precursor mass tolerance, while the CID data used 0.6 m/z fragment tolerance and the ETD data used 1.2 m/z fragment tolerance. All searches used fixed carbamidomethylation, variable phosphorylation (S,T,Y), oxidation (M), and pyro-Glu (peptide N-term Q). Finally, the ptmRS node was appended to the end of the workflow to calculate modification site localization probability.

The Consensus workflow is used to calculate protein groups, filtering peptides and proteins by false discovery rate, and rolling up the quantification results from individual peptide spectral matches (PSM's) to peptide groups and proteins.



FIGURE 3. Study factors for the enriched (a) phosphopeptide search and the flow-through and wash fractions (b). There are as many columns in the final report as are shown in green on this table.

The final step is to set up the columns to be displayed in the final result for quantification. The columns that are highlighted in green as in Figure 3 are those that will show up in the final report with the quantification values. This setup also works similarly for reporter ion quantification and isotopically-labeled precursor quantification.

For these data, two different analyses were run. The first analyzed the samples from the enriched fractions to compare phosphopeptides while the second analyzed the combined flow-through and wash samples to compare changes in protein abundances across the samples.

Results

Phosphopeptide search results

The IMAC enrichment was less successful for these samples and thus the IMAC enrichment data for samples 3955 and 3956 were removed from the analysis for clarity. As a result, there were two technical replicates for the TiO_2 -enriched phosphopeptides for each of the original samples leading to 8 different quantitative categories as shown in Figure 4 for the 242 raw files. Figure 4 shows a screenshot of the Proteome Discoverer results of the phosphopeptide data.



Figure 4. Identified peptide groups for TiO₂-enriched phosphopeptides. The fourth column shows the modification and the site localization probability calculated by ptmRS. The "Area columns" show the peak area calculated for that identified phosphopeptide across all 8 multidimensional LC/MS/MS runs. The selected peptide is about 20x more abundant in the hESCs compared to the hNSCs with good biological and technical reproducibility.

Table 1 shows a summary of the number of PSM's, peptides, and phosphopeptides identified across the 4 samples. To find phosphorylation sites that are unique to the two hESC samples (3955 and 3957), the row filters were selected to show phosphopeptides that appear in one of the two replicates in each case that were not identified in the two hNSC samples (3956 and 3958). In total, there were 811 unique phosphopeptides identified only in the hESC samples versus 253 unique phosphopeptides identified only in the two hNSC samples.

Sample	PSM's	Unique peptides	Unique phoshpopeptides	Unique phosphopeptides (ptmRS isoform score >60)						
hESC 1	38174	7715	5575	4151						
hNSC 1	21987	4562	3191	2388						
hESC 2	47338	7835	6223	4526						
hNSC 2	35350	7086	4898	3677						
TABLE 1 Identification summary for enriched phosphopentide samples										

TABLE 1. Identification summary for enriched phosphopeptide samples.

For the unique hESC phosphopeptides, overrepresentation analysis of pathways was performed and insulin signaling, MAPK, ErbB, and AMPK pathways were identified. For the peptides unique to the hNSC samples, there were several phosphorylation sites detected on the protein MAP2, which is known to be involved in neurogenesis, SOX-5, which is involved in chondrogenesis, and neuron-navigator 1. Some phosphopeptides were also identified in all samples but are also differentially abundant. There were 38 phosphopeptides that appear to be up-regulated by a factor of ≥ 2 in hNSCs while there were 101 phosphopeptides that were up-regulated in hESCs. The phosphopeptides with the highest differential expression ratios between the hESC and hNSC samples are shown in Table 2.

Peptide	Protein	Ratio hNSC/hESC
LKCGSGPVHISGQHLVAVEEDAE <u>S</u> EDEEEEDVK	Rho GTPase activating protein 17	0.025
RPTPNDDTLDEGVGLVHSNIATEHIP S PAK	Hydroxymethylglutaryl-CoA synthase	0.032
MAPTPIPTR S PSDSSTASTPVAEQIER	Drebrin	0.055
SSM <u>S</u>GLHLVK	Acetyl-CoA carboxylase 1	0.058
DME <u>S</u> PTKLDVTLAK	Microtubule associated protein 4	0.068
TTR <u>T</u> PEEGGYSYDISEK	Microtubule associated protein 1B	18
RPA <u>S</u> PSSPEHLPATPAESPAQR	Sin3 histone deacetylase corepressor complex component SDS3	19
VAL <u>S</u> DDETKETENMR	DNA polymerase delta subunit 3	21
RS <u>T</u> QGVTLTDLQEAEK	Protein phosphatase 1 regulatory subunit 12A	31
IED <u>S</u> EPHIPLIDDTDAEDDAPTKR	Plasma membrane calcium-transporting ATPase 1	120

TABLE 2. Phosphopeptides with the lowest and highest quantitative ratios between hESCs and hNSCs.

Protein Quantification

For protein quantification, another set of searches were performed on the flow-through and wash fractions from the phosphopeptide enrichment steps. The same workflows were used as shown in Figure 2, but phosphorylation was removed as a variable modification from the Sequest HT searches and Protein FDR threshold node set to 1% was added. For these data, the "Proteins" tab in the final results show the average of the top 3 most abundant peptides detected across all of the samples as seen in Figure 5 below.

g.	Check	ed Maste	er Accession	Description	Coverage	f Papidas	s PSHe	# Unique Pepfides	8 Protein Groupe	1 AA	M-/ JiCu	calc pl	2005. T.C.C. Plan decorphismed. 1	2055, T.C.C. Plan Broughinson, 2	2005, ISBN: Flue Besudylinesh, 2 2005, ISBN: Flue Besudylinesh, 2	2006, ISBC: Plan Encodylensis, 1	2006, 19IAC, Plan Ewoldhissoh, 2	2006, T.C.C. Flow throughingsh. 1	2006, 1022, First Broughtman, 2 2007, TDD, First Broughtman, 1	2807, TGC, Plan Broughinstei, 2	2008, TCO, Flow Broughtmach, 1	2008, T.CO. Pars Recognised 2	Found is	in Sa
0		1	P42704	Laucine rich FPR motif containing protein, mitochondrial	65%	25	2229	25	1	1354	157.8	6.13	9247	9547	24 13	61 9247	9747	9247 1	Ce3 1174	0 1940	1041	15+5		
		1Ý	075083	htD repeat containing protein 1	77%	45	1798	45	1	606	88.2	6.65	1.845	1.845	545 2.74	1 324	3.445	2745 2	045 3.5e	8 3745	2445	2545	27	11
5 4		1V	P22234	multifunctional protein ADE2	623	43	2103	43	1	425	47.0	7.23	2.245	2245 0	5el 30e	4 33el	3.745	6345 E	245 4 14	4.424	4.642	4.845	27	11
5 -0		1¥	P15311	Earin	69%	54	2348	29	1	500	69.4	6.27	1545	1540	Sel 204	4 12.7	8.047	1241 1	40 24	4 2848	1248	1248	17	11
7 0		V	Q57457-2	Isoform 2 of E3 ubiquitir-protein ligase UBR4	275	140	2165	145	1	5204	\$75.6	6.01	2347	3547 4	447 5.24	3 4567	5047	5047 5	047 5.04	0 6247	6167	6.6e7		11
1.11		V	OFFICE 1	Hypoxia up-regulated protein 1	561	87	2054	32	1	999	111.3	5.22	4.347	6.847 0	447 5.24	Titel	1741	I fel 1	141 1 1 1	5 11al	2148	2345		
		12	Q14315	Filamin-C	525	114	2091	22	1	2725	250.8	5.97	7247	8 147 4	547 5.04	0 5947	6 147	9.347 9	67 8 14	0 840	5747	5 447	27	11
		1 Ý	Q12931	heat shock protein 75 kDa. mitochondrial	551	47	2117	48	1	704	80.1	8.21	1.5+5	1341	343 14	0 2140	2349	1542 1	10 24	0 2540	1040	2040	27	11
1.0		1V	P46777	60S ribosomal protein L5	62%	30	1952	33	1	297	34.3	9.72	3.5+5	35-8	044	4 514	5.5+5	62+5 6	545 5.74	4 5646	6748	57+5	27	11
		V	Q9P258	Protein ROC2	225	35	2044	36	1	522	56.0	8.78	1345	1340	1ad 3.94	4 3.6+5	33+5	2848 2	Nel 3.04	4 2940	2.040	2.348		11
1		1 V	QRUDRS	Alcha-aninoadipic semisidehude sunthase mitochondrial	58%	63	2177	61	1	125	102.1	6.64	9.547	1141.00	fel 164	1.1.1	144	1241 1	al 15	1.174	15al	1 fel		
		V	P17907	T-complex protein 1 subunit siphs	685	41	2164	41	1	556	60.3	6.11	1.64	1041	Pel 214	4 204	234	4341 4	141 H 44	1.534	3.64	274		
5 44		14	P62325	OTP-binding nuclear protein RAN	563	25	2003	25	1	216	24.4	7.49	4.045	4 245	Cel 10	1445	1549	5.145 5.		8 8745	1245	1249		
		1Ý	Q99832	T-complex protein 1 subunit ets	71%	48	2292	19	1	543	55.3	7.65	1.745	1240 11	541 2.34	6 2146	2240	35+5 2	940 2.24	0 2040	2440	22+0	27	11
7 -		V	P06753-2	Isoform 2 of Tropomyosin alpha-3 chain	62%	35	1788	3	1	248	29.0	4.78	8.447	8.547	1647 5.04	0 4547	4.747	7.847 7	547 8.54	0 940	8.167	8.047		11
1. 10		1¥	P27816-1	Microtubule associated protein 4	66%	75	2160	31	1	1152	120.9	5.43	4.547	5.447 1	447 504	0.000	1.545	1545 1	145 144	4 204	8.547	0.5e7	17	11
		V	P49508	Atanine-6R50A ligase, cytoplasmic	8122	56	1935	56	1	260	106.7	5.53	9247	9347	Sel 16	4 124	1248	1241 1	348 2.24	4 224	124	124		11
5 -0		V	Q01225	transpelin	54%	28	2117	28	1	201	22.6	0.04	4.540	4340 0	442 5.94	4 224	2440	2340 2		6 1045	2540	2748		11
		V	P23395-1	405 ribosomal protein 53	89%	40	2260	40	1	243	267	2.00	5240	5.540 0	345 5.64	4 7.045	7.545	0.140 0.	140 11.54	6 116	3040	9245		6.8
			Q13283	Res GTPase activating protein-binding protein 1	675	27	1809	28	1	466	52.1	5.52	1545	1545	548 24	4 15e5	1.545	2545 2	45 2.54	8 2.545	2345	2245		
		1V	P30041	Perovinedoxin-6	73%	25	2062	26	1	224	25.0	6.18	3.045	3248 3	1al 35	4.444	\$ 245	8345 8	148 2.14	4 144	1.545	8.348		
		1¥	09+972	Protein lin-28 homolog A	785	1 13	1963	19	1	229	22.7	8.07	2345	2545 3	18el 38e	4 134	7.565	9.445 9		4 424	1248	1 245	17	11
5 0		V	P07237	Protein disutide-isomerase	69%	43	1971	43	1	500	57.1	4.07	1240	1240 11	Sel 15	4 294	32+0	2548 2	040 174	6 1740	2140	2748		11
		V	P11505	C-1-tetrahydrofolate synthese, cytoplasmic	65%	67	1961	66	1	905	101.5	7.30	1.440	1540	548 254	1.645	1040	1540 1	548 2.84	2 2040	1540	1545		11
7 -14		V	P63244	Ousnine nucleotide-binding protein subunit beta-2-like 1	240	33	1996	33	1	317	35.1	7.65	3.645	3745 1	345 5.04	4 1041	6.445	6345 5	545 S 14	5 5040	5440	5 545	6 R	
		ΤÝ	P07226-1	Tropomyosin alpha-4 chain	82%	27	1681	20	1	248	28.5	4.69	9747	104	245 1.24	4 144	1.445	2.445 2	145 1.54	0.176	2140	2245		1
		Y	Q15233	Non-POU domain-containing octamer binding protain	68%	43	2069	40	1	471	54.2	8.95	1545	1741.0	1al 24	4 354	1141	5248 5	64 23	0 2040	1141	140	17	0
5 0		1¥	P63104-1	14-3-3 protein zetaldelta	845	30	2160	23	1	245	27.7	4.79	5240	5240 1	741 5.74	2445	9.440	1.50 1	00 110	0 110	1049	1049		11
1 -		V	Q954E4	Far upstream element-binding protein 1	68%	43	2122	3	1	644	67.5	7.61	2.547	2247 4	1547 454	4767	\$ 147	2447 2	347 3.24	0 120	4947	4.547		11
2		V	Q15851	UTP-glucose-1-phosphate uridylyfosnelerase	77%	50	2154	50	1	506	56.2	8.15	2540	3148 4	045 4 14	4 254	4.545	2545 3		6 524	1040	3748		11
3. 44		V	Q14340-2	Isoform 2 of Eukaryotic initiation factor 43-8	88%	31	2043	13	1	406	46.5	5.48	4.547	4.547	1947 4.34	d 11el	1.145	1045 1	145 7.84	0 7.507	1140	1245	6 R	6.8
4		ΤÝ	P05023	Sodumbotassium-transporting ATPase subunit alpha-1	525	54	1801	32	1	1023	112.8	5.49	1.1+0	1240 0	041 204	6 7.8+7	8547	1540 1	40 22	0 2240	1540	1540	- 7	11

Ready 10062 Protein Group: 10052/24319 Protein: 10050 Peptide Group:: 1704732 PSMI: 305165 MS-MS Spectrum Into: 125 Result

Figure 5. Identified proteins for the flow-through and wash fractions across the samples. The "Areas" table shows the average of the peak areas of the top 3 identified peptides from that protein for each sample.

A total of 10682 unique proteins were identified across all of the samples. Table 3 shows the proteins identified for each of the samples, with roughly 6500-7000 proteins identified in each sample. Row filters were applied to show only those proteins identified in the various hESC samples, with a total of 285 proteins with two or more unique peptides identified in at least one of the hESC samples but none of the hNSC samples (data not shown). This includes proteins such as cadherin-3, PR domain zinc finger protein 14, Tyrosine-protein kinase Lck, and Oct4. A second set of filters were used to show proteins that only appear in one or more hNSC samples, which produced a list of 458 proteins ephrin type-A receptor 4 isoform a precursor, contactin-associated protein 1 precursor, and Spondin-1.

Sample	Enrichment	Replicate	Protein groups	Peptide Groups
Sample	Enrichment	Replicate	Protein groups	Peptide Groups

hESC 1	TiO ₂	1	6240	47667
hESC 1	TiO ₂	2	6303	48105
hESC 1	IMAC	1	5854	35924
hESC 1	IMAC	2	5991	36990
hNSC 1	TiO ₂	1	7047	54935
hNSC 1	TiO ₂	2	7044	55473
hNSC 1	IMAC	1	6551	43585
hNSC 1	IMAC	2	6654	44475
hESC 2	TiO ₂	1	6701	54023
hESC 2	TiO ₂	2	6776	54385
hNSC 2	TiO ₂	1	6605	50303
hNSC 2	TiO ₂	2	6639	50887

TABLE 3. Summary of protein identifications for each sample.

The table of protein identifications and quantitative information was exported to Excel. To account for differences in the injection amounts for each sample, the abundances for each protein were normalized to the summed abundance of all proteins in the sample. The results were imported into ProteinCenter and profiled. The top and bottom clusters showed overrepresentation in hNSCs and hESCs respectively (Figure 8).



FIGURE 8. Top two clusters from ProteinCenter profiling. The upper cluster corresponds to proteins more abundant in the hNSCs whereas the lower cluster corresponds to proteins that are more abundant in the hESCs.

Some selected proteins in cluster 1 include several ephrin receptors, PAX-6, and frizzled-3 precursor all of which are involved in neuronal development in animal models, and SOX-2, which is known to be involved in stem cell pluripotency and differentiation. Cluster 2 contains dozens of highly up-regulated (≥10x) proteins including gamma-synuclein, CD9 antigen, calveolin-1 isoform alpha, HRAS-like suppressor 3, and nocturnin.

Conclusion

- Proteome Discoverer 2.0 software is well equipped to analyze highly complex datasets, in this case a dataset with well over 800 RAW data files.
- The study management feature was used to produce peak area quantification values for phosphopeptides and the proteins from the flow-through and wash fractions from TiO₂ and IMAC enrichment.
- The precursor ion quantification can be used to find differentially expressed phosphopeptides as well as proteins between the hESC and hNSC samples.
- ProteinCenter aids in the biological interpretation of the Proteome Discoverer software results.

www.thermoscientific.com

©2015 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Denmark +45 70 23 62 60 Europe-Other +43 1 333 50 34 0 Finland +358 10 3292 200 France +33 1 60 92 48 00 Germany +49 6103 408 1014 India +91 22 6742 9494 Italy +39 02 950 591 Japan +81 45 453 9100 Korea +82 2 3420 8600 Latin America +1 561 688 8700 Middle East +43 1 333 50 34 0 Netherlands +31 76 579 55 55 New Zealand +64 9 980 6700 Norway +46 8 556 468 00



 Russia/CIS
 +43 1 333 50 34 0

 Singapore
 +65 6289 1190

 Spain
 +34 914 845 965

 Sweden
 +46 8 556 468 00

 Switzerland
 +41 161 716 77 00

 UK
 +44 1442 233555

 USA
 +1 800 532 4752

