

# Standards Chemical Kit

With Low/High Concentration PPGs (PN 4406127)



Insert PN 4412466 Rev. B  
Printed in USA

For Research Use Only.  
Not for use in diagnostic procedures.

For all chemicals, read the MSDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## List of kit components

Item (ingredients)	Volume (mL)	Quantity (each)	Applied Biosystems part number
POS PPG, <sup>†</sup> 1E-4 M (100 µM or stock solution) [Polypropylene glycol, average M <sub>n</sub> ~425 <sup>‡</sup> ; polypropylene glycol, average M <sub>n</sub> ~1000; polypropylene glycol, average M <sub>n</sub> ~2000; ammonium acetate; formic acid; ACS reagent, 96%; acetonitrile, HPLC-grade; methanol; HPLC-grade; water]	20	1	4405227
POS PPG, 1E-5 M (10 µM or 10:1) <sup>§</sup>	20	1	4405229
POS PPG, 2E-6 M (2 µM or 50:1) <sup>§</sup>	20	1	4405231
POS PPG, 2E-7 M (0.2 µM or 500:1) <sup>§</sup>	20	1	4405233
NEG PPG, 3E-4M (300 µM or stock solution) [Polypropylene glycol, average M <sub>n</sub> ~425; polypropylene glycol, average M <sub>n</sub> ~1000; polypropylene glycol, average M <sub>n</sub> ~2000; polypropylene glycol, average M <sub>n</sub> ~3000; ammonium acetate; formic acid; ACS reagent, 96%; acetonitrile; HPLC-grade; methanol; HPLC-grade; water]	20	1	4405234
NEG PPG, 3E-5 M (30 µM or 10:1) <sup>††</sup>	20	1	4405235
RESERPINE, 1 pmol/µL (1 µM) [Reserpine, acetonitrile, acetic acid, water]	5	1	4405236
RESERPINE, 0.167 pmol/µL (0.167 µM or 6:1) <sup>††</sup>	5	2	4405237
RESERPINE, 0.0167 pmol/µL (0.0167 µM or 60:1) <sup>††</sup>	5	2	4405238
ES TUNING MIX <sup>§§</sup> [Betaine; trifluoroacetic acid ammonium salt; hexamethoxyphosphazene; hexakis(2,2-difluoroethoxy)Phosphazene; hexakis(1H, 1H, 3H-tetrafluoropropoxy)phosphazene; hexakis(1H, 1H, 5H-octafluoropentoxy)phosphazene; hexakis(1H, 1H, 7H-dodecafluoroheptoxy)phosphazene; hexakis(1H, 1H, 9H-perfluorononyloxy)phosphazene; tris (heptafluoropropyl)-S-triazine; acetonitrile; distilled-deionized water]	4	1	4405243
AMMONIUM ACETATE [Ammonium acetate; renin substrate tetradecapeptide; porcine; acetic acid, HPLC-grade]	2 g	1	4405245
Renin-Reserpine Mix (1 pmol/µL Renin + 400 fmol/µL Reserpine) [Reserpine; Acetonitrile, HPLC-grade; renin substrate tetradecapeptide; porcine; acetic acid; water]	5	1	4405247
WASH SOLVENT [50/50 methanol/water]	30	1	4405249
REN/RES DILUTION SOLVENT [50/50 acetonitrile/water + 0.1% acetic acid]	20	1	4405255
ES TUNING DILUTION SOLVENT [95/5 acetonitrile/water]	20	1	4412397
BOTTLE, HDPE, CLEAR NM RND, 15 mL	—	1	T8183

<sup>†</sup> Poly(propylene glycol)

<sup>‡</sup> Narrow molecular-weight distribution

<sup>§</sup> Same ingredients as in POS PPG, 1E-4 M (100 µM)

<sup>††</sup> Same ingredients as in NEG PPG, 3E-4 M (300 µM)

## Same ingredients as in RESERPINE, 1 pmol/μL (1 μM)

§§ Prepare as described in "100:1 ES Tuning Mix."

When stored at 2 to 8 °C, this kit maintains performance for one year. Before use, mix components well.

## Prepare solutions

### 100:1 ES Tuning Mix

Mix in the provided empty bottle 9.9 mL of ES TUNING DILUTION SOLVENT and 100 μL ES TUNING MIX stock solution. The diluted solution can be used for up to one week after preparation if the diluted solution is stored at 2 to 8 °C.

### Reconstituted stock solutions

If any stock solution requires reconstitution, it is recommended to use the reconstituted solution within one week after preparation while the solution is stored at 2 to 8 °C.

### Solvents

Prepare additional solvents, if required:

Preparation of solvents†

Component	POS PPG Dilution Solvent	NEG PPG Dilution Solvent	REN/RES DILUTION SOLVENT	WASH SOLVENT	ES TUNING DILUTION SOLVENT	Flow Injection Solvent 70/30 ACN/water‡
Methanol	50 mL	75 mL	—	50 mL	—	—
Formic acid	100 μL	100 μL	—	—	—	0.1 mL
Acetonitrile	100 μL	100 μL	50 mL	—	95 mL	700 mL
Acetic acid	—	—	0.1 mL	—	—	—
Ammonium acetate	15.4 mg	391 mg	—	—	—	154 mg
Water	49.8 mL	24.8 mL	49.9 mL	50 mL	5 mL	299.9 mL

† All the chemicals for preparing dilution solvents should be HPLC-grade or higher.

‡ Required for selected instruments only.

Store the diluted solvent at 2 to 8 °C.

## How to obtain support

For the latest services and support information for all locations, go to:

[www.appliedbiosystems.com](http://www.appliedbiosystems.com), then click **Support**.

At the Support home page, you can:

- Access worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.
- Submit a question directly to Technical Support.
- Order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents.

© Copyright 2009, Applied Biosystems. All rights reserved.

**For Research Use Only. Not for use in diagnostic procedures.**

NOTICE TO PURCHASER: LIMITED LICENSE

Applied Biosystems, and AB (Design) are Registered trademarks of Applied Biosystems or its subsidiaries in the US and/or certain other countries.

All other trademarks are the sole property of their respective owners.