

Standards Chemical Kit With Low/High Concentration PPGs

Product No. **4406127**
Lot No. **0910050**
Date of Manufacture **02 Nov 2009**
Expiration Date: **10 Jun 2010**

Concentration			
BOTTLE PART NUMBER	BULK PART NUMBER	COMPONENT	CONCENTRATION (+/- ≤2%)
4405227	4405198	POS PPG 1E-4M	100 µM
4405229	4406096	POS PPG 1E-5M	10 µM
4405231	4406098	POS PPG 2E-6M	2 µM
4405233	4406100	POS PPG 2E-7M	0.2 µM
4405234	4405200	NEG PPG 3E-4M	300 µM
4405235	4406102	NEG PPG 3E-5M	30 µM
4405236	4405202	RESERPINE 1 µM	1 µM
4405237	4406104	RESERPINE 0.167 µM	0.167 µM
4405238	4406106	RESERPINE 0.0167 µM	0.0167 µM
4405247	4405208	RENIN-RESERPINE MIX	1 µM - 0.4 µM
(1 pmol/µL Renin + 400 fmol/µL Reserpine)			

Assurance of Purity and Traceability

These standards have been manufactured using ISO9001 procedures. Concentrations are calculated gravimetrically, and balances are calibrated to NIST traceable standards in accordance with our ISO procedures. Purity from interferences is verified through mass spectrometric analysis. Solvents are minimum HPLC grade.

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

**ISO9001
REGISTERED**

Foster City, CA
Pleasanton, CA
Bedford, MA
Austin, TX
Warrington, UK
Singapore



Quality Assurance Manager

To obtain a Certificate of Analysis on-line go to www.appliedbiosystems.com or email us at cofarequest@appliedbiosystems.com