



## Product Information Sheet

Product Name: Peptide Mass Standard Kit

Catalog Number: AS-60882

Size: 1Kit

Description: This kit consists of two calibration mixtures for calibrating mass scale in MALDI-TOF or ESI mass spectrometry. Calibration Mixture 1 ranges from 800 to 1800 daltons, and Calibration Mixture 2 from 1800 to 3800 daltons.

### Standard Operation Protocol:

1. Reconstitute Calibration Mixture 1 and Calibration Mixture 2 each in 100  $\mu$ l of Standard Diluent (Component C).
2. Prepare matrix solution by dissolving 1 vial of Component D in 100  $\mu$ l of matrix diluent (Component E).
3. Mix 24  $\mu$ l of matrix solution with 1  $\mu$ l of Calibration Mixture 1 or 1  $\mu$ l of Calibration Mixture 2 to obtain Matrix: Standard solutions. Load 1  $\mu$ l per sample position.

### Kit Components:

Component A: Calibration Mixture 1 (1 vial)

Component B: Calibration Mixture 2 (1 vial)

Component C: Standard Diluent (30% Acetonitrile in 0.1% TFA, 250  $\mu$ l)

Component D: Matrix ( $\alpha$ -Cyano-4-hydroxycinnamic acid, 2 vials)

Component E: Matrix Diluent (50% Acetonitrile in 0.1% TFA, 1 ml)

### Calibration Mixture 1:

Name	(M+H) <sup>+</sup> Average	(M+H) <sup>+</sup> Monoisotopic	Size
Des-Arg1-Bradykinin	905.05	904.47	10 $\mu$ g, (+/-)10%
Angiotensin I	1297.51	1296.68	40 $\mu$ g, (+/-)10%
Neurotensin	1673.96	1672.92	10 $\mu$ g, (+/-)10%

### Calibration Mixture 2:

Name	(M+H) <sup>+</sup> Average	(M+H) <sup>+</sup> Monoisotopic	Size
ACTH (Clip 1-17)	2094.46	2093.09	10 $\mu$ g, (+/-)10%
ACTH (Clip 18-39)	2466.72	2465.2	40 $\mu$ g, (+/-)10%
ACTH (Clip 7-38)	3660.19	3657.923	40 $\mu$ g, (+/-)10%



**Storage:** Store all kit components at -20°C. Long term storage of reconstituted peptide is not recommended.

Please note that the peptide solution is not stable. We recommend using it within 24-48 hours of dissolution.