

Batch information

STD

HTRF Hu. Amyloid b1-42 kit - Std

Product reference :

62B42CDA**Product characteristics :**

| | |
|-----------------------|---|
| - Lot Number | 01A |
| - Conditioning buffer | Phosphate 50 mM pH 7.0 + BSA 1% + stabilizers |
| - Appearance | Lyophilized form |
| - Storage conditions | 2-8 °C |
| - Expiration date | 2024.02.22 |

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (MAb) | 1 800.00 pg | N/A |

PLEASE NOTE

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

