

Batch information

STD HTRF Human IL4 kit - Std

Product reference : **62HIL04CDA**

Product characteristics :

| | |
|-----------------------|--|
| - Lot Number | 02A |
| - Conditioning buffer | Phosphate 50 mM pH7.0 + BSA 1% + stabilizers |
| - Appearance | Lyophilized form |
| - Storage conditions | 2-8 °C |
| - Expiration date | 2024.12.05 |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 4 200.00 pg | N/A |

PLEASE NOTE

Reconstitute with 0.350 mL H₂O.

The expiration date of batch 02A has been extended to 2024-12-05 after validation by QC.

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

