



Batch information

STD

HTRF Human IL23 kit - Std

Product reference :

62HIL23CDA

Product characteristics :

- Lot Number 04A
- Conditioning buffer Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance Lyophilized form
- Storage conditions 2-8 °C
- Expiration date 2025.08.17

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	31 500.00 pg	N/A

PLEASE NOTE

Reconstitute with 0.350 mL H2O.

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