

STD

Human IL12 kit - Standard

Product reference :

62HIL12CDA**Product characteristics :**

- Lot Number_____ 04B
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 1% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.09.11

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	15 600.00 µg	N/A

PLEASE NOTEReconstitute with 0.400 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

