

Human IL-11 Standard

Human IL11 kit - Standard

Product reference :

62HIL11CDA

Product characteristics :

- Lot Number_____ 01A
- Conditioning buffer_____ Phosphate 50mM pH7.0 +BSA 1% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.11.25
- Active Moiety Concentration (STD)_____ 15000 pg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (STD)	3 150.00 pg	N/A

PLEASE NOTE

Reconstitute with 210µl 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

