

Batch information

Human IL-11 Standard

HTRF Human IL11 kit - Std

Product reference :

62HIL11CDA**Product characteristics :**

- Lot Number	02A
- Conditioning buffer	Phosphate 50mM pH7.0 +BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.10.27
- Active Moiety Concentration (STD)	15000.00 pg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (STD)	4 500.00 pg	N/A

PLEASE NOTEReconstitute with 300µl 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

