



Technical Data Sheet

STD

HTRF Mouse IL6 kit - Std

Product reference :

62MIL06CDA**Product characteristics :**

- Lot Number	05AR
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2026.11.07

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	6 300.00 pg	N/A

PLEASE NOTE

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