

Batch information

cIAP1 BIR2 Standard

HTRF cIAP1 BIR2 Bind Kit - Std

Product reference :

64BDIAPB2CDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	100% DMSO
- Appearance	Frozen form
- Storage conditions	-16 °C
- Expiration date	2024.06.22
- Volume per vial	30.00 µL
- Active Moiety Concentration (STD)	20.00 mM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (STD)	600.00 nmoles	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

