

Batch information

cIAP1 BIR3 Standard

HTRF cIAP1 BIR3 Bind Kit - Std

Product reference :

64BDIAPB3CDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	Tris Hepes 50 mM pH7 + 0.1 % BSA + stabilizer
- Appearance	Frozen form
- Storage conditions	-16 °C
- Expiration date	2024.06.22
- Volume per vial	30.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (std)	3.00 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

