

Batch information

IBMX

HTRF IBMX - 40 µL

Product reference :

62AMXADA

Product characteristics :

- Lot Number	05A
- Conditioning buffer	DMSO
- Appearance	Liquid form
- Storage conditions	4 °C
- Expiration date	2024.11.25
- Active Moiety Concentration (0)	500.00 mM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (0)	40.00 µl	222

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

