

Batch information

Control lysate

HTRF GAPDH House. kit - Ctrl lysate

Product reference :

64GAPDHTDA

Product characteristics :

- Lot Number	04A
- Conditioning buffer	Lysis Buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2025.02.03
- Volume per vial	150.00 µL
- Active Moiety Concentration (CL)	1.00 X

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (CL)	150.00 µl	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

