

Batch information

Control lysate HTRF H2AX p-S139 kit - Ctrl lysate
Product reference : **64H2XTDA**

Product characteristics :

- Lot Number 11A
- Conditioning buffer Cellular kinase lysis buffer
- Appearance Frozen form
- Storage conditions -60 °C
- Expiration date 2025.03.07
- Volume per vial 150.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	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

