

## Batch information

**VHL binding kit std**

HTRF Human VHL Bind. Kit - Std

Product reference :

**64BDVHLCDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	HEPES 50 mM pH7.0 + BSA 0.1% + stabilizers
- Appearance	Frozen form
- Storage conditions	-16 °C
- Expiration date	2024.12.01

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (STD)	50.00 nmoles	473

### 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](http://www.cisbio.com)*

