

## Batch information

**LCL161 Red Ligand**

HTRF LCL161 Red Ligand

Product reference :

**64BDIAPB3RED**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	Tris Hepes 50 mM Ph7+ 0.1% BSA + 0.05% Tween20 + Stabilizer
- Appearance	Frozen form
- Storage conditions	-16 °C
- Expiration date	2024.06.22
- Volume per vial	200.00 µL
- Active Moiety Concentration (Lig)	8.00 µM

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Lig)	1.60 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](http://www.cisbio.com)*

