

**D2V2-NHS**

d2v2-NHS - 1 mg

Product reference :

**65DVSABB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 07A  
- Conditioning buffer \_\_\_\_\_ ///  
- Appearance \_\_\_\_\_ Desiccated form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2023.04.20

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (DV)	1.00 mg	1 066

**PLEASE NOTE**

Rz ( D0650nm/280nm)=34.7

Reconstitute with DMSO

***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)***