

## CERTIFICATE OF ANALYSIS

**Catalog Number** C049  
**Product Name** NBD C<sub>6</sub>-ceramide,  
[6-((N-(7-Nitrobenz-2-oxa-1,3-diazol-4-yl)amino)caproyl) sphingosine]  
**CAS Number** 86701-10-2  
**Molecular Formula** C<sub>30</sub>H<sub>49</sub>N<sub>5</sub>O<sub>6</sub>  
**Molecular Weight** 575.75  
**Appearance** orange solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	466 nm 23,600 cm <sup>-1</sup> M <sup>-1</sup>	466 ± 4 nm 22,500 ± 2,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup> Emission Maximum	537 nm	534 ± 4 nm
<b>TLC</b> Purity	96 %	≥ 90 %
<b>NMR</b> Result	meets specification	consistent with structure

1. Solvent: Methanol.