

NBD C6-Sphingomyelin

NBD C6-sphingomyelin (6-((N-(7-nitrobenz-2-oxa-1,3-diazol-4-yl)amino)hexanoyl)sphingosyl phosphocholine) has been used for following sphingolipid metabolism in cells.



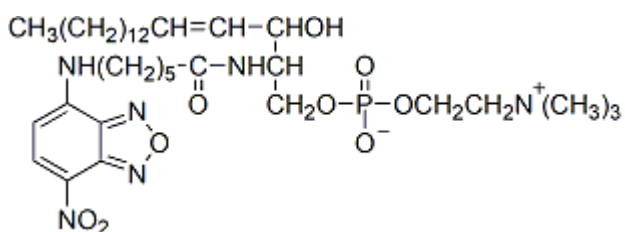
Product attributes

CAS number	94885-04-8
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Green
Excitation/Emission	466/536 nm
Conjugation	NBD

Product Description

NBD C6-sphingomyelin (6-((N-(7-nitrobenz-2-oxa-1,3-diazol-4-yl)amino)hexanoyl)sphingosyl phosphocholine) has been used for following sphingolipid metabolism in cells.

- $\lambda_{\text{Ex}}/\lambda_{\text{Em}}$ (MeOH) = 466/536 nm
- Dark solid soluble in chloroform or DMSO
- Store at -20 °C and protect from light
- $\text{C}_{35}\text{H}_{61}\text{N}_6\text{O}_9\text{P}$
- MW: 740.87
- [94885-04-8]



References

1. Methods Cell Biol. 38, 221 (1993).
2. Biochim Biophys Acta 1113, 277 (1992).
3. Adv Cell Mol Biol Membranes 1, 199 (1993).
4. Biochim Biophys Acta 1082, 113 (1991).

This datasheet was generated on June 12, 2026 at 02:01:25 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nbd-c6-sphingomyelin-6-n-7-nitrobenz-2-oxa-1-3-diazol-4-ylamino-hexanoyl-sphingosyl-phosphocholine/>