



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Pluronic® F-127

Nonionic detergent useful for solubilizing hydrophobic compounds to facilitate cell loading.



Product attributes

CAS number	9003-11-6
------------	-----------

Product Description

Pluronic® F-127 is a nonionic detergent useful for solubilizing hydrophobic dye molecules such as cell-permeable AM esters of Fura-2, Fluo-3, Indo-1, or Rhod-2 calcium indicators to facilitate cell loading. Pluronic® F-127 is usually dissolved in water or DMSO with gentle heating. We also offer ready-to-use solutions of [Pluronic® F-127, 20% in DMSO](#) and [Pluronic® F-127, 10% in H₂O](#).

- White solid soluble in water and DMSO
- Store at room temperature
- MW: ~12,500
- [9003-11-6]

Pluronic F-127 is a registered trade mark of BASF.

References

- 1) J. Membrane Biol. 19, 1 (1974).
- 2) Science 233, 886 (1986).
- 3) J. Biol Chem 262, 12801 (1987).
- 4) J. Biol. Chem. 265, 19543 (1990).
- 5) Methods Enzymol. 302, 341 (1999).
- 6) Methods Enzymol. 307, 441(1999).

This datasheet was generated on January 15, 2026 at 12:27:42 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/pluronic-f-127/>