

Pluronic® F-127, 10% in H2O

Ready-to-use solution of Pluronic® F-127, a nonionic detergent useful for solubilizing hydrophobic compounds like AM esters of calcium indicator dyes to facilitate cell loading.



Product attributes

Call us: 800-304-5357

CAS number

9003-11-6

Product Description

Ready-to-use solution of Pluronic® F-127, a nonionic detergent useful for solubilizing hydrophobic dye molecules such as AM esters of Fura-2, Fluo-3, Indo-1, or Rhod-2 to facilitate cell loading.

We also offer Pluronic® F-127 solid and Pluronic® F-127, 20% Solution in DMSO.

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

References

- 1) J. Membrane Biol. 19, 1 (1974).

- 1) J. Melhibaria 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 November 26, 2025 at 04:48:35 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/pluronic-f-127-10-in-h2o/