

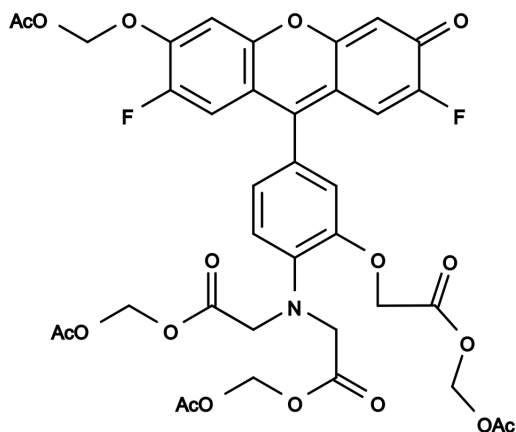
## Product Information Sheet

### Ordering Information

Product Number:	20401
Product Name:	Mag-Fluo-4 AM
Unit Size:	10x50 ug
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	817.65
Chemical Structure:	



Soluble In:	DMSO
Excitation Wavelength:	494
Emission Wavelength:	516

### Application Notes

The cell-permeant Mag-Fluo-4 AM is an analog of Fluo-4 AM with a  $K_d$  for Mg ion of 4.7 mM and a  $K_d$  for Ca ion of 22  $\mu$ M, making it useful as an intracellular Mg ion indicator as well as a low-affinity Ca ion indicator. This low-affinity fluorescent Ca ion indicator has been used to accurately track the kinetics of the spatially averaged free Ca ion transient in skeletal muscle. Mag-fluo-4 yields reliable kinetic information about the spatially averaged free Ca ion transient in skeletal muscle.