Fluo-3 AM (ultra pure)

Calcium dye

Fluo-3 AM is widely used in flow cytometry, confocal laserscanning microscopy and fluorescence spectroscopy. The probe has proved particularly valuable for determining GPCR function via cell-based high-throughput screening. Fluo-3 AM is a membrane permeable probe and can be passively loaded in cells by simple incubation. Once inside the cell, Fluo-3 becomes fluorescent in the presence of free Ca (II) . Fluo-3 is a long wavelength calcium indicator probe (488 nm excitable). The longer wavelength of absorption is convenient for minimizing photodamage in cells. Wavelength Maxima: Excitation 506nm, Emission 526nm

Citations: 4

View Online »

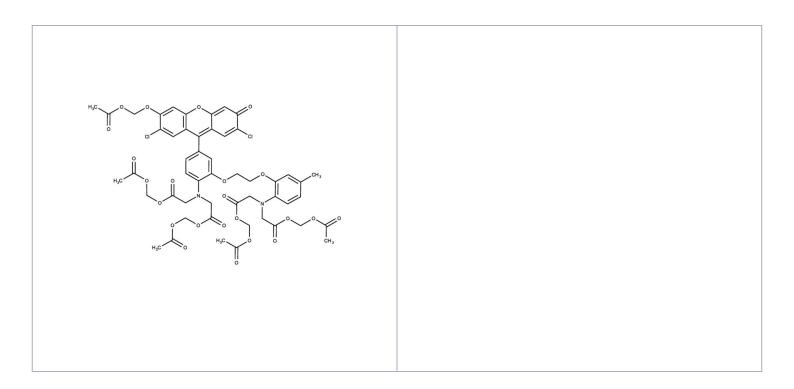
Ordering Information

Order Online »

ENZ-52004 1mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability Stable for at least one year after receipt when stored as recommended.

Handling Protect from light. Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-[Amino-5-(2,7-dichloro-6-acetomethoxy-3-oxo-3H-

xanthen-9-yl)phenoxy]-2-(2'-amino-5'-

methylphenoxy)ethane-N,N,N',N'-tetraacetic acid,

pentaacetoxy-methyl ester

CAS 121714-22-5

MW 1129.8

Purity ≥95% (HPLC)

Solubility Soluble in DMSO.

Technical Info / Product NotesThis product is a member of the CELLESTIAL[®] product

line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL® reagents and kits are optimal for use in

demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency

and reproducibility are required.

Last modified: May 29, 2024

