

# PRODUCT INFORMATION



## Fluo-3 (sodium salt)

Item No. 20404

**Formal Name:** N-[2-[2-[2-[bis(carboxymethyl)amino]-5-(2,7-dichloro-6-oxido-3-oxo-3H-xanthen-9-yl)phenoxy]ethoxy]-4-methylphenyl]-N-(azanediyl)-diacetate, pentasodium salt

**MF:** C<sub>36</sub>H<sub>25</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>13</sub> • 5Na

**FW:** 879.4

**Purity:** ≥90%

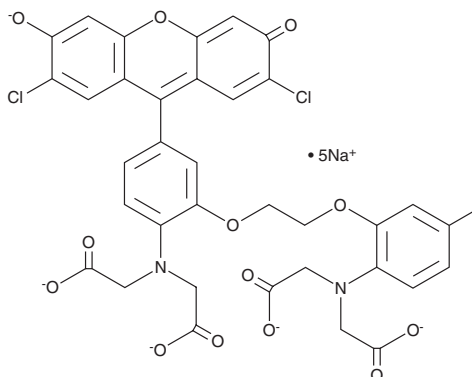
**UV/Vis.:** λ<sub>max</sub>: XX nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥2 years

**Special Conditions:** Protect from light



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Fluo-3 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the fluo-3 (sodium salt) in the solvent of choice. Fluo-3 (sodium salt) is soluble in water.

### Description

Fluo-3 is a fluorescent calcium indicator commonly used in flow cytometry and cell-based experiments to detect changes in intracellular calcium levels.<sup>1</sup> Its absorption maximum at 506 nm makes it compatible with excitation at 488 nm by argon-ion laser sources. Fluo-3 provides intense fluorescence upon binding calcium, detected at a maximum emission at 526 nm which can be monitored by FL1 (green, 525 nm band pass) sensors in flow cytometry.

### Reference

1. Minta, A., Kao, J. P., and Tsien, R. Y. Fluorescent indicators for cytosolic calcium based on rhodamine and fluorescein chromophores. *J. Biol. Chem.* **264**(14), 8171-8178 (1989).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/11/2021

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM