

# PRODUCT INFORMATION



## Indo-1 (sodium salt)

Item No. 20419

**Formal Name:** 1-[2-amino-5-(6-carboxyindol-2-yl)phenoxy]-2-(2'-amino-5'-methylphenoxy)ethane]-N,N,N',N'-tetraacetic acid, pentasodium salt

**MF:** C<sub>32</sub>H<sub>26</sub>N<sub>3</sub>O<sub>12</sub> • 5Na

**FW:** 759.5

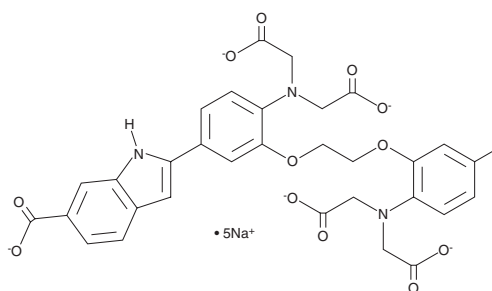
**Purity:** ≥90%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Special Conditions:** Light sensitive



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

### Laboratory Procedures

Indo-1 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the indo-1 (sodium salt) in the solvent of choice, which should be purged with an inert gas. Indo-1 (sodium salt) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Indo-1 is a ratiometric fluorescent calcium indicator. It is ideal for analyses using flow cytometry, as it uses a single excitation source, typically 349-364 nm light from an argon-ion laser. The emission maximum shifts from 475-485 nm without calcium to 400-410 nm when indo-1 binds calcium.<sup>1,2</sup> Indo-1 is prone to photobleaching, which limits its usefulness in methods involving microscopy.<sup>1</sup>

### References

1. Paredes, R. M., Etzler, J. C., Watts, L. T., *et al.* Chemical calcium indicators. *Methods* **46(3)**, 143-151 (2008).
2. Grynkiewicz, G., Poenie, M., and Tsien, R. Y. A new generation of Ca<sup>2+</sup> indicators with greatly improved fluorescence properties. *J. Biol. Chem.* **260(6)**, 3440-3450 (1985).

#### 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, 10/17/2022

#### 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