

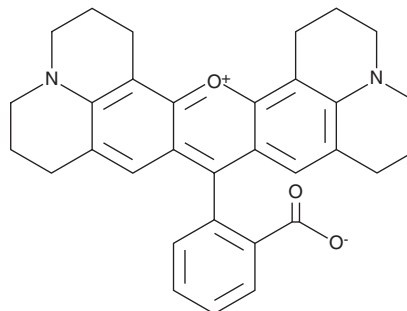
# PRODUCT INFORMATION



## Rhodamine 101 (inner salt)

Item No. 20654

**CAS Registry No.:** 116450-56-7  
**Formal Name:** 9-(2-carboxyphenyl)-2,3,6,7,12,13,16,17-octahydro-1H,5H,11H,15H-xantheno[2,3,4ij:5,6,7-i'j'] diquinolizin-18-ium, inner salt  
**Synonyms:** RH101, Rhodamine 640  
**MF:** C<sub>32</sub>H<sub>30</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 490.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 232, 266, 311, 367, 567 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Rhodamine 101 (inner salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the rhodamine 101 (inner salt) in the solvent of choice, which should be purged with an inert gas. Rhodamine 101 (inner salt) is soluble in the organic solvent ethanol at a concentration of approximately 0.10 mg/ml.

### Description

Rhodamine 101 (inner salt) is a bright fluorescent dye with excitation and emission maxima of 565 and 595 nm, respectively.<sup>1,2</sup> It can be used in various biological applications such as fluorescence microscopy, flow cytometry, fluorescence correlation spectroscopy, and ELISA.

### References

1. Aigner, D., *et al.* New fluorescent perylene bisimide indicators--a platform for broadband pH optodes. *Anal. Bioanal. Chem.* **400(8)**, 2475-2485 (2011).
2. Crissman, H. A., and Steinkamp, J. A. Rapid, one step staining procedures for analysis of cellular DNA and protein by single and dual laser flow cytometry. *Cytometry* **3(2)**, 84-90 (1982).

#### 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, 12/08/2022

#### 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