

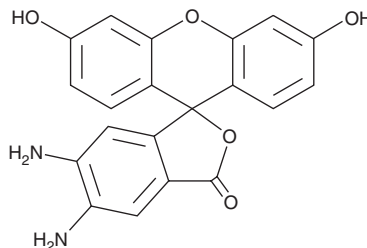
PRODUCT INFORMATION



DAF-2

Item No. 85160

CAS Registry No.: 205391-01-1
Formal Name: 2-(3,6-dihydroxy-4,5-diamino-9H-xanthen-9-yl)-benzoic acid
Synonyms: 4,5-Diaminofluorescein
MF: C₂₀H₁₄N₂O₅
FW: 362.3
Purity: ≥98%
Ex./Em. Max: 485/538 nm
Supplied as: A solution in DMSO
Storage: -20°C
Stability: ≥2 years



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

Description

DAF-2 is a sensitive fluorescent indicator commonly used for the detection of nitric oxide (NO).^{1,2} It reacts with NO in the presence of oxygen to yield the highly fluorescent triazolofluorescein (DAF-2T). Fluorescence is monitored using excitation and emission wavelengths of 485 and 538 nm, respectively.² At neutral pH the detection limit for NO is 2-5 nM.

References

1. Kojima, H., Sakurai, K., Kikuchi, K., *et al.* Development of a fluorescent indicator for nitric oxide based on the fluorescein chromophore. *Chem. Pharm. Bull. (Tokyo)* **46(2)**, 373-375 (1998).
2. Nakatsubo, N., Kojima, H., Kikuchi, K., *et al.* Direct evidence of nitric oxide production from bovine aortic endothelial cells using new fluorescence indicators: Diaminofluoresceins. *FEBS Lett.* **427(2)**, 263-266 (1998).

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, 12/13/2022

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