

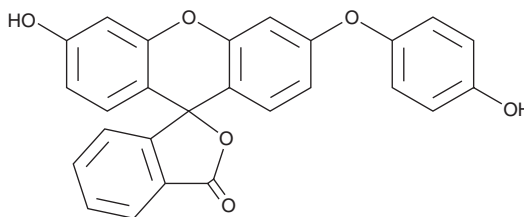
PRODUCT INFORMATION



HPF

Item No. 10159

CAS Registry No.: 359010-69-8
Formal Name: (2-[6-(4'-hydroxy)phenoxy-3H-xanthene-3-on-9-yl]benzoic acid
MF: C₂₆H₁₆O₆
FW: 424.4
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 277 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

HPF is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the HPF under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of HPF in these solvents is approximately 20 mg/ml.

HPF is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of HPF should be diluted with the aqueous buffer of choice. HPF has a solubility of approximately 0.15 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method.

Description

HPF is a cell-permeable aromatic amino-fluorescein derivative that has little intrinsic fluorescence.¹ It undergoes oxidation only by highly reactive oxygen species (hROS) such as the hydroxyl radical, peroxyxynitrite, and hROS generated from a peroxidase/H₂O₂ system. It is inert to hypochlorite ion, nitric oxide, hydrogen peroxide (H₂O₂), superoxide, and other oxidants. Upon oxidation, HPF is converted to the highly fluorescent molecule fluorescein, with excitation/emission maxima of 490/515 nm, respectively, allowing the simple direct detection of highly reactive biological radicals.

Reference

1. Setsukinai, K.i., Urano, Y., Kakinuma, K., *et al.* Development of novel fluorescence probes that can reliably detect reactive oxygen species and distinguish specific species. *J. Biol. Chem.* **278**(5), 3170-3175 (2003).

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, 02/01/2019

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