

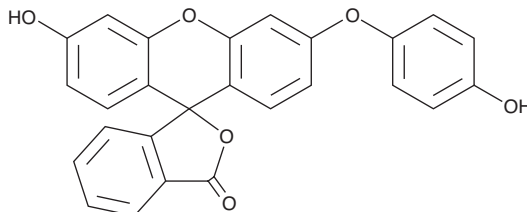
Product Information



HPF

Catalog No. 10159

Formal Name: (2-[6-(4'-hydroxy)phenoxy-3H-xanthene-3-on-9-yl]benzoic acid
MF: C₂₆H₁₆O₆
FW: 424.4
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that HPF be stored as supplied at -20°C. It will be stable for at least one year.

HPF is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate 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 methyl acetate 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. We do not recommend storing the aqueous solution for more than one day.

The biology of highly reactive oxygen radical species is of great interest in many biomedical research disciplines, including neurodegeneration and aging, cancer, and infectious diseases.¹ There are a number of fluorescent reagents, such as 2,7-dichlorodihydro-fluorescein (DCDHF), that can be used to detect free radicals, but they have significant limitations due to their facile oxidation by light and numerous non-radical oxidants such as hydrogen peroxide (H₂O₂).² HPF is a cell-permeable aromatic amino-fluorescein derivative that has little intrinsic fluorescence. It undergoes oxidation only by highly reactive oxygen species (hROS) such as the hydroxyl radical, peroxynitrite, and hROS generated from a peroxidase/H₂O₂ system. It is inert to hypochlorite ion, nitric oxide, H₂O₂, superoxide, and other oxidants. Upon oxidation, HPF is converted to the highly fluorescent molecule fluorescein, allowing the simple direct detection of highly reactive biological radicals.³

References

1. Matés, J.M., Pérez-Gómez, C., and Nuñez de Castro, I. Antioxidant enzymes and human diseases. *Clin. Biochem.* **32**(8), 595-603 (1999).
2. Hempel, S.L., Buettner, G.R., O'Malley, Y.Q., *et al.* Dihydrofluorescein diacetate is superior for detecting intracellular oxidants: Comparison with 2',7'-dichlorodihydrofluorescein diacetate, 5-(and 6)-carboxy-2',7'-dichlorodihydrofluorescein diacetate, and dihydrorhodamine 123 *Free Radic. Biol. Med.* **27**(1), 146-159 (1999).
3. Setskunai, K., 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).

Related Products

APF - Cat. No. 10157 • DAN-1 EE (hydrochloride) - Cat. No. 85070 • Dihydrorhodamine 123 - Cat. No. 85100 • 2,7-Dichlorodihydrofluorescein diacetate - Cat. No. 85155 • DAF-2 diacetate - Cat. No. 85165

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 03/21/2006