

APF

Fluorescent probe for hROS detection

Fluorescent reagent (Ex(max): 490nm; Em(max): 515nm) for the detection of highly reactive oxygen species (hROS). Immediately reacts with hROS such as hydroxyl radical, peroxynitrite and hypochlorite, and the fluorescence intensity greatly increases. Use of APF together with HPF (Prod. No. ALX-620-074) also allows for specific detection of hypochlorite (OCl^-) to elucidate reliable the roles of OCl^- in biological systems such as neutrophils. In addition, peroxynitrite can be detected in distinction from nitric oxide and superoxide since APF does not react with nitric oxide, superoxide and hydrogen peroxide. Moreover, APF is resistant to light-induced autooxidation. *Not for sale in Japan.*

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-620-075-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

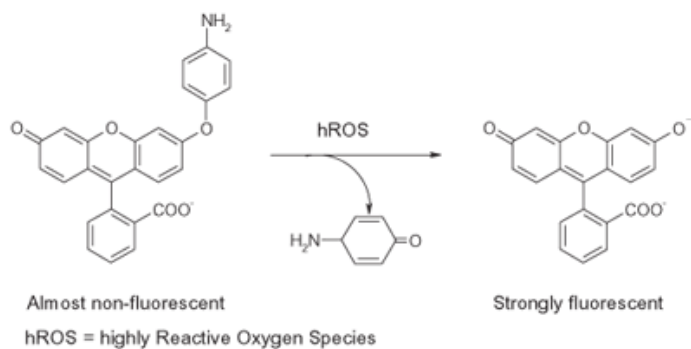


Figure 1: Reaction of Aminophenyl Fluorescein (APF) with hROS.

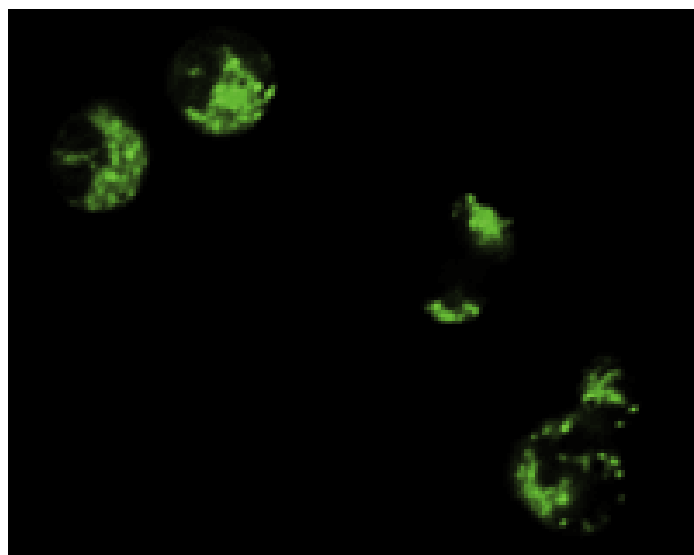
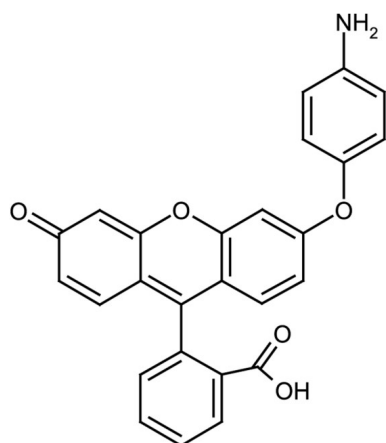


Figure 6: Fluorescence image of hypochlorite (OCl^-) which was generated in porcine neutrophils using APF.



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare 500-5'000-fold dilution (~10-1µM) in phosphate buffer (0.1M phosphate, pH 7.4) immediately before use. BSA, phenol red and amines may affect the fluorescence and must be used with caution. Do not store the dilutions.
Handling	After opening, prepare aliquots and store at +4°C. Protect from light. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Aminophenyl fluorescein, 2-[6-(4'-Amino)phenoxy-3H-xanthen-3-on-9-yl]benzoic acid
Appearance	Pale yellow liquid.
CAS	359010-70-1
Formula	$C_{26}H_{17}NO_5$
Formulation	Dissolved in 0.47ml dimethyl formamide.
MW	423.4
Purity	≥98% (HPLC)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)