

5-Aminosalicylic acid

NF-κB pathway inhibitor

Inhibits TNF-α regulated IκB degradation and NF-κB activation.

Citations: 2

[View Online »](#)

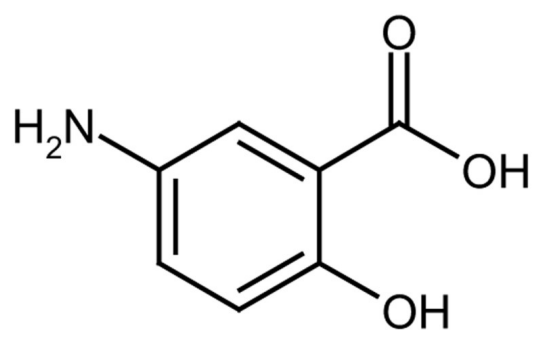
Ordering Information

[Order Online »](#)

ALX-430-110-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	5-ASA, Mesalamine, 5-Amino-2-hydroxybenzoic acid
Appearance	White to light gray or light pink powder or crystalline solid.
CAS	89-57-6
Formula	$C_7H_7NO_3$
Identity	Determined by IR.
MI	14: 5904
MW	153.1
Purity	≥98% (TLC)
Solubility	Soluble in 1 M hydrochloric acid.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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