

5-AIQ . hydrochloride

PARP-1 inhibitor

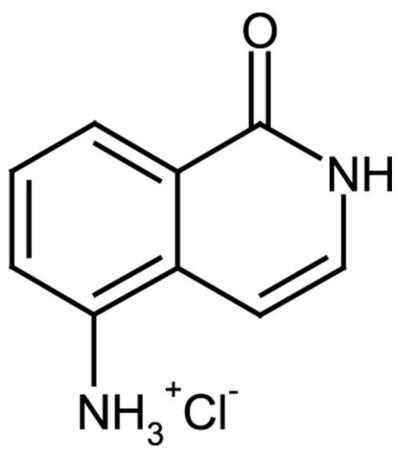
Water-soluble, potent inhibitor of poly(ADP-ribose) polymerase-1 (PARP-1).

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Aminoisoquinolinone . HCl
Appearance	Beige to brown powder.
CAS	93117-07-8
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_9H_8N_2O$. HCl
MW	160.2 . 36.5
Purity	≥97% (HPLC)
Solubility	Soluble in water (warm).

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