

Quinolinic Acid monoclonal antibody

Ordering Information

[Order Online »](#)

IMS-IS002	50µl
-----------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store at +4°C for short term (1-2 months). Aliquot and store at -20°C for long term. Avoid repeated freeze / thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Pyridine-2,3-dicarboxylic acid, 2,3-pyridinedicarboxylic acid, 3,4-Pyridinedicarboxylic acid, Pyridine-3,4-dicarboxylic acid
Application	IF, IHC
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Conjugated Quinolinic acid
Isotype	IgG1 k chain
Recommendation Dilutions/Conditions	Immunohistochemistry (IHC) – Dilute at 1:100-1:1000 Immunofluorescence (IF) – Dilute at 1:50-1:500
Species Reactivity	Species independent
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



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