

Protein Phosphatase 2A1 (bovine kidney), (purified)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	PP2A1
Application Notes	Useful for enzyme kinetics studies, drug screening.
Formulation	Liquid. In 50 mM Tris, pH 7.0, containing 0.1 mM EDTA, 50% glycerol, 14 mM 2-mercaptoethanol, 1 mM benzamidine, 0.1 mM PMSF.
MW	36 55 60 kDa
Source	Isolated from bovine kidney.

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