

ADP-ribosylation factor monoclonal antibody (1D9)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ARF

Application Flow Cytometry, ICC, IF, IHC, WB

Application Notes Detects a band of ~21kDa by Western blot.
Not suitable for neutralizing ARF in solution.

Clone 1D9

Formulation Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Recombinant human ARF1 (ADP-ribosylation factor-1).

Isotype IgG1

Recommendation Dilutions/Conditions Flow Cytometry (1:20)Immunocytochemistry / Immunofluorescence (1:100)Immunohistochemistry (1:100)Western Blot (1:3000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Dog, Hamster, Human, Mouse, Rat

Specificity Recognizes ARF1, ARF3, ARF5 and ARF6.

UniProt ID P84077

Last modified: May 29, 2024



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)