

ADP-ribosylation factor monoclonal antibody (1D9)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-028-R100	100µl
------------------	-------

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 Flow Cytometry (1:20)Immunocytochemistry / Immunofluorescence
Dilutions/Conditions (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



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-be@enzolifesciences.com

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

France
Phone: +33 1 72 44 01 555
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