

NF-ATc1 monoclonal antibody (7A6)

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

Citations: 27

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-0220	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 NFATc1, Nuclear factor of activated T cells, cytoplasmic 1, NFAT2, NFATC

Application CHIP, Gel Shift, ICC, IF, IHC, IP, WB

Application Notes Detects a band of ~105kDa by Western blot.

Clone 7A6

Crossreactivity Does not cross-react with NF-ATc2.

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

Host Mouse

Immunogen Bacterially expressed glutathione S-transferase (GST) fusion protein containing NF-ATc1 residues 1-654.

Isotype IgG1

Recommendation Dilutions/Conditions Immunocytochemistry (1:250)Immunofluorescence (1:10-1:100)Immunohistochemistry (1:20-200)Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Monkey, Mouse, Rat

Technical Info / Product Notes

Note: Confusion may arise regarding NFAT protein versus gene names. According to UniProt, the nomenclature is as follows:**Nuclear Factor of Activated T Cells, Cytoplasmic 1 (NF-ATc1)** Gene Names: NFATC1 or NFATC or NFAT2

Nuclear Factor of Activated T Cells, Cytoplasmic 2 (NF-ATc2) Gene Names: NFATC2 or NFATP or NFAT1

Nuclear Factor of Activated T Cells, Cytoplasmic 3 (NF-ATc3) Gene Names: NFATC3 or NFATX or NFAT4

Nuclear Factor of Activated T Cells, Cytoplasmic 4 (NF-ATc4) Gene Names: NFATC4 or NFAT3

Nuclear Factor of Activated T Cells (NF-AT5) Gene Names: NFAT5

UniProt ID

O95644

Worry-free Guarantee

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

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