

NF-ATc2 monoclonal antibody (25A10.D6.D2)

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-215-C100	100µg
------------------	-------

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 NFATc2, Nuclear factor of activated T cells, cytoplasmic 2, NFATP, NFAT1

Application Gel Supershift, ICC, IF, IHC (FS), IHC (PS), IP, WB

Application Notes Detects bands of ~140kDa (phosphorylated NF-ATc2 in resting immune cells) and ~120kDa (fully-dephosphorylated NF-ATc2 in stimulated cells) by Western blot.

Clone 25A10.D6.D2

Crossreactivity Does not cross-react with NF-ATc1.

Formulation Liquid. Purified antibody in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 51-69 (A51ISSPSGLAYPDDVLDYGL69) of mouse nuclear factor of activated T cells, cytoplasmic 2 (NF-ATc2) conserved in all three isoforms A, B and C.

Isotype IgG1

Recommendation Dilutions/Conditions Immunocytochemistry (1:100 – 1:1,000)Immunofluorescence (1:100 – 1:1,000)Immunohistochemistry (frozen sections and paraffin sections, 1:50 – 1:200)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, 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 Q60591

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



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)