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.

lsotype lgG1

Recommendation Dilutions/Conditions Immunocytochemistry (1:100 –

1:1,000)lmmunofluorescence (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.

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

Worry-free Guarantee

Last modified: May 29, 2024

Q60591

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, Phone: 800.942.0430 infousa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com

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

Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900

uk@enzolifesciences.com