

# NF-ATc1 monoclonal antibody (7A6)

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

Citations: 27

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-022-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	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](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)