

Tollip monoclonal antibody (Kimmy-2)

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

Ordering Information

[Order Online »](#)

ALX-804-154-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

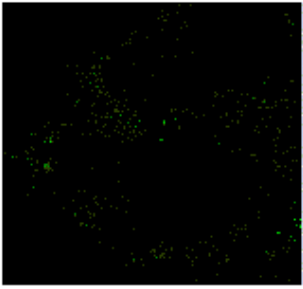
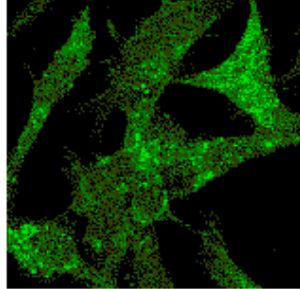
A**B**

Figure: Immunocytochemistry showing the detection of Tollip in Tollip -/- (A) or overexpressed Tollip in -/- (B) murine embryonic fibroblasts (MEF) cells with the MAb

to Tollip (Kimmy-2) (Prod. No. [ALX-804-154](#)).**Method:** Following fixation with paraformaldehyde 3.7% and permeabilization with 0.2% saponin, MEF cells were

incubated with the MAb to Tollip (Kimmy-2) at a 1:100 dilution. An anti-mouse IgG antibody conjugated with Alexa Fluor 488 was used for detection.”””,Tollip monoclonal antibody (Kimmy-2)

Immunocytochemistry,Tollip monoclonal antibody (Kimmy-2) Immunocytochemistry,yes

ALX-804-157,804-

157.gif,[http://static.enzolifesciences.com/fileadmin/files/image/804-](http://static.enzolifesciences.com/fileadmin/files/image/804-157.gif)

157.gif,**Method:** HEK 293T cells were mock transfected (thin line) or transfected with an expression plasmid enabling surface expression of the extracellular domain of mouse BAFF (thick line). Cells (5×10^5) were incubated on ice for 30 min in 50 μ l FACS buffer (PBS

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ICC, IP, WB
Clone	Kimmy-2
Crossreactivity	Weakly cross-reacts with human tollip.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant mouse Tollip (aa 1-274).
Isotype	IgG1
Recommendation Dilutions/Conditions	Immunocytochemistry (1:100)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue supernatant.
Species Reactivity	Mouse
UniProt ID	Q9QZ06
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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