

BTLA monoclonal antibody (BTLA/1528)

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

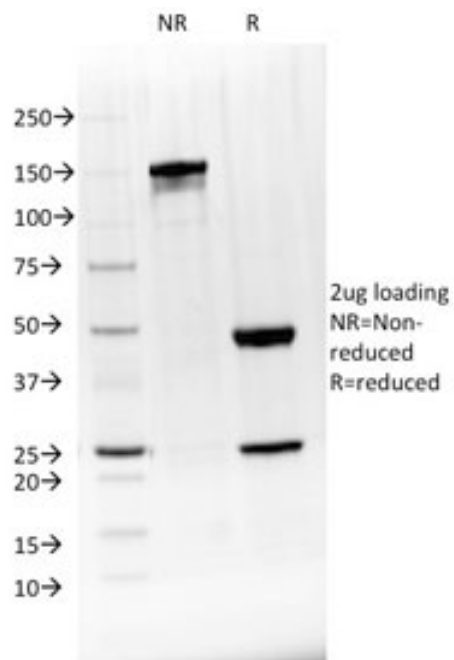
Ordering Information

[Order Online »](#)

ENZ-ABS871-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of purified, BSA-free BTLA antibody (clone BTLA/1528) as confirmation of integrity and purity.

Handling & Storage

Handling	Store the BTLA antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide).
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	B- and T-lymphocyte attenuator, B- and T-lymphocyte-associated protein, CD272
Application	ELISA
Clone	BTLA/1528
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant extracellular human CD272/BTLA.
Isotype	IgG1κ
Purity Detail	Protein G affinity chromatography purified.
Recommendation Dilutions/Conditions	Optimal conditions must be determined individually for each application.
Species Reactivity	Human
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-be@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)