

# BTLA monoclonal antibody (BTLA/1528)

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

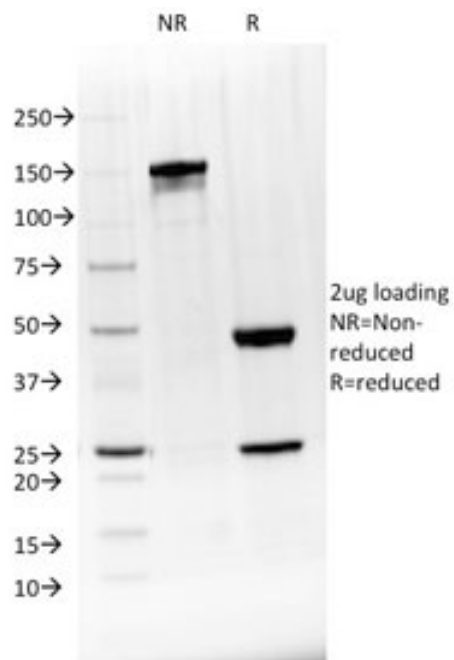
## Ordering Information

[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 <a href="#">Worry-Free Guarantee</a>