

ST2 polyclonal antibody

ST2 is a member of a superfamily containing the interleukin-1 receptor and the Toll-like receptors (TLRs). ST2 was originally identified as a protein whose production was stimulated by various proliferation inducing agents such as PDGF and FGF. Additionally, it has been shown to negatively regulate IL-1 receptor and Toll-like receptor (TLR) 4 signaling and to maintain endotoxin tolerance.

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

Citations: 1

[View Online »](#)

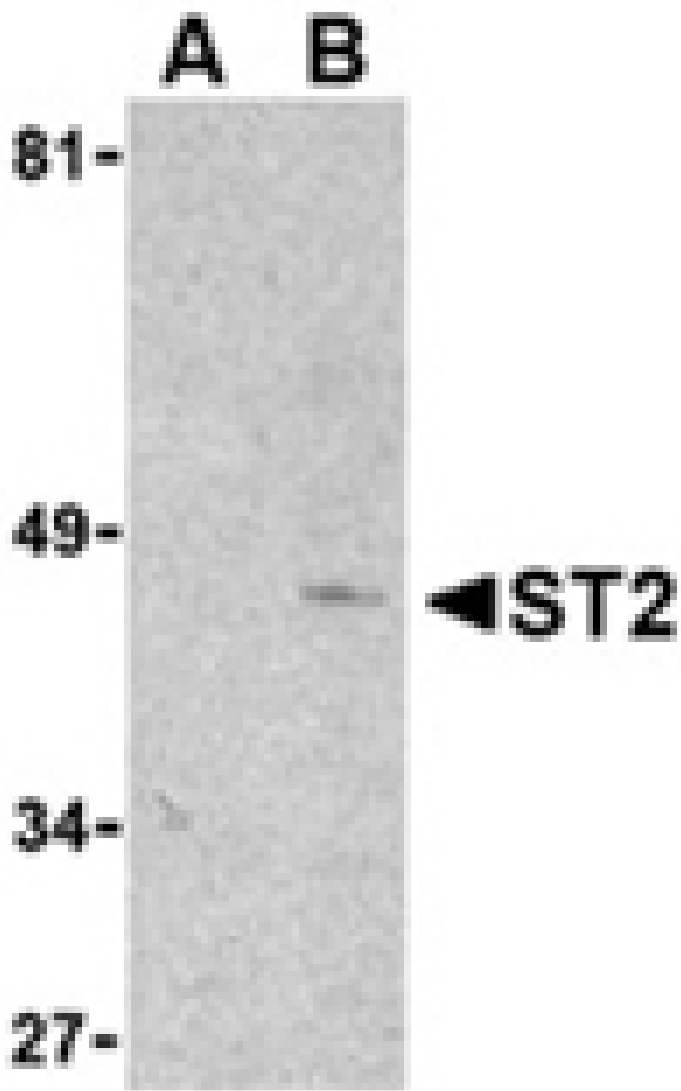
Ordering Information

[Order Online »](#)

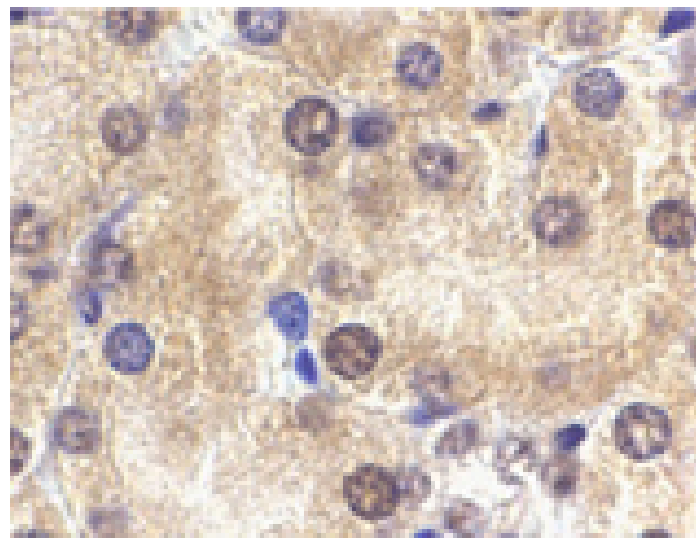
| | |
|---------------|-------|
| ADI-CSA-512-E | 100µg |
|---------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of ST2 in mouse kidney lysate with ST2 antibody at 1 μ g/ml in the presence (lane A) or absence (lane B) of 1 μ g blocking peptide.



Immunohistochemistry analysis of ST2 in mouse kidney tissue with ST2 antibody at 2 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------------------------|---|
| Alternative Name | Interleukin-1 receptor-like 1, Lymphocyte antigen 84 |
| Application | IHC, WB |
| Application Notes | Detects a band of ~45kDa by Western blot. |
| Formulation | Liquid. In PBS containing 0.02% sodium azide. |
| GenBank ID | BAA02854 |
| Host | Rabbit |
| Immunogen | Synthetic peptide corresponding to the sequence near the N-terminus of mouse ST2. The sequence is common to all three ST2 isoforms. |
| Purity Detail | Peptide affinity purified. |
| Recommendation Dilutions/Conditions | Immunohistochemistry (2µg/ml)Western Blot (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity | Human, Mouse |
| UniProt ID | P14719 |
| Worry-free Guarantee | This antibody is covered by our Worry-Free Guarantee |



ENZO LIFE SCIENCES,
INC.
Phone: +1 940 948 9120
info@enzolifesciences.com
usa@enzolifesciences.com

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

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

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

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

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com