

Anti-dsRNA monoclonal antibody K1 (500 µg)

SCICONS Anti-dsRNA monoclonal antibody K1 can be used for the detection of dsRNA both in cultured cells and fixed paraffin-embedded histological samples. It is recommended for Poly I:C detection (higher affinity than J2) and SCICONS Anti-dsRNA monoclonal antibody J2 alternative to resolve cross-reactions and/or remove background in those rare experimental setups where J2 does not provides satisfactory results. Applications: ELISA, IF, FACS, IHC, IP, Dot Blot, ChIP, affinity purification, immunoelectron microscopy. Please note that nucleic acid separation prior to dsRNA-immunoblotting must be carried out by polyacrylamide gel electrophoresis, because the sensitivity of detection is considerably lower after blotting from agarose gels.

Ordering Information

[Order Online »](#)

JBS-RNT-SCI-10020500	500µg
----------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation lyophilized from a 1 mg/ml solution in PBS

Host Mouse

Isotype IgG2a, kappa

Reconstitution To prepare a 1 µg/µl PBS antibody solution add 200 µl (RNT-SCI-10020200) or 500 µl (RNT-SCI-10020500) sterile DNase/RNase-free distilled water. As a result of the lyophilisation procedure, the reconstituted antibody may contain small amounts of denatured protein in the form of aggregates that may interfere with some applications such as immunohistochemistry (e.g. by giving high backgrounds). We therefore highly recommend to spin down (microcentrifuge) the reconstituted antibody before use and to use the supernatant only.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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)