

# Anti-dsRNA monoclonal antibody K2 (10 ml)

SCICONS Anti-dsRNA monoclonal antibody K2 can be used for the detection of dsRNA. It is recommended for (Sandwich-) ELISA and provides an isotype alternative to SCICONS Anti-dsRNA monoclonal antibodies J2 and K1. Applications: (Sandwich-)ELISA (1:2 dilution in PBS) , IHC, Dot Blot

## Ordering Information

[Order Online »](#)

JBS-RNT-SCI-10030010	10ml
----------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Formulation** undiluted hybridoma cell culture supernatant (RPMI 1640)

**Host** Mouse

**Isotype** IgG2a, kappa

**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)