



Anti-dsRNA monoclonal antibody K2

mouse, IgM, kappa chain

Cat. No.	Amount
RNT-SCI-10030005	5 ml
RNT-SCI-10030010	10 ml



For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C to -80 °C upon reconstitution for long-term storage

Additional Storage Conditions: avoid freeze/thaw cycles, store in aliquot.

Shelf Life: 12 months after date of delivery

Form: undiluted hybridoma cell culture supernatant (RPMI 1640)

Description:

Anti-dsRNA monoclonal antibody K2 can be used for the detection of dsRNA. It is recommended for (Sandwich-) ELISA and provides an isotope alternative to Anti-dsRNA monoclonal antibodies J2 and K1.

Applications:

(Sandwich-)ELISA (1:2 dilution in PBS) , IHC, Dot Blot

Specificity:

Anti-dsRNA monoclonal antibody K2 recognises double-stranded RNA (dsRNA) provided that the length of the helix is greater than or equal to 40 bp dsRNA. Recognition is independent of the sequence and nucleotide composition of the antigen. All naturally occurring dsRNAs investigated up to now (40-50 species) as well as poly(I)-poly(C) and poly(A)-poly(U) have been recognised by Anti-dsRNA monoclonal antibody K2.

As described by Schönborn et al. (1991), Anti-dsRNA monoclonal antibody K2 binds with high avidity to all dsRNAs investigated.

Species Origin: Mouse

Heavy Chain Isotype: IgM

Light Chain Isotype: kappa

Quality control:

Activity: AN-ELISA (relative activity compared to reference K2)

Selected References:

Schönborn et al. (1991) Monoclonal antibodies to double-stranded RNA as probes of RNA structure in crude nucleic acid extracts. *Nucleic Acids Res.* **19**: 2993.

Lukacs (1994) Detection of virus infection in plants and differentiation between coexisting viruses by monoclonal antibodies to double-stranded RNA. *J. Virol. Methods* **47**: 255.

Lukacs (1997) Detection of sense:antisense duplexes by structure-specific anti-RNA antibodies. In: *Antisense Technology. A Practical Approach*, C. Lichtenstein and W. Nellen (eds), pp. 281-295. IRL Press, Oxford