

# dsRNA Recombinant Monoclonal Mouse Antibody (rJ2), Lyophilized - Biotium Choice



Recombinant J2 antibody for detection of double-stranded RNA (dsRNA). Suitable for IHC, IF, ELISA, EMSA, flow cytometry, and dot blot.

## Product Description

dsRNA Recombinant Monoclonal Mouse Antibody (rJ2) detects double-stranded RNA (dsRNA). This antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance. This antibody is suitable for use in IHC, IF, ELISA, EMSA, flow cytometry, and dot blot. Each lot is tested to verify that it is RNase-free.

dsRNA Recombinant Monoclonal Mouse Antibody (rJ2) recognizes double-stranded RNA (dsRNA) as long as the helix is at least 40 base pairs in length. Antigen recognition is independent of the RNA sequence or nucleotide composition. The J2 antibody has been validated by numerous publications and is shown to detect naturally occurring dsRNAs, including synthetic dsRNAs such as poly(I)-poly(C) and poly(A)-poly(U). However, in some assays, its affinity for poly(I)-poly(C) is approximately 10-fold lower than for other dsRNA antigens.

**Reconstitution:** Add nuclease-free dH<sub>2</sub>O to the vial to dissolve the antibody at 1 mg/mL (200 uL for 200 ug vial or 500 uL for 500 ug vial). Mix gently to dissolve completely. After reconstitution, antibody concentration will be 1 mg/mL in 1X PBS with no BSA or azide.

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise for flow cytometry and other applications. [View our full catalog of Biotium Choice antibodies.](#)

## References

1. Nucleic Acids Res. 1991 Jun 11;19(11):2993-3000. DOI: [10.1093/nar/19.11.2993](https://doi.org/10.1093/nar/19.11.2993).

This datasheet was generated on June 15, 2026 at 02:04:12 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/dsrna-recombinant-monoclonal-mouse-antibody-rj2-lyophilized-biotium-choice/>

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

<b>Antibody number</b>	#P025
<b>Biotium Choice Antibody</b>	Primary
<b>Antibody type</b>	Biotium Choice, Primary
<b>Clonality</b>	Monoclonal, Recombinant Monoclonal
<b>Host species</b>	Mouse
<b>Clone</b>	rJ2
<b>Isotype</b>	IgG2a, kappa
<b>Antibody reactivity (target)</b>	dsRNA
<b>Synonyms</b>	Double-stranded RNA; dsRNA
<b>Antibody target cellular localization</b>	Cytoplasm
<b>Verified antibody applications</b>	IF (verified), Dot blot (verified)
<b>Positive control</b>	poly(I-C) dot blot, HeLa cells
<b>Antibody application notes</b>	Stains endogenous dsRNA in HeLa cells by immunofluorescence. Selectively detects dsRNA on dot blot.
<b>Antibody research areas</b>	RNA Analysis & Detection, Vaccine Research, Immunology, Virology
<b>Antibody/conjugate formulation</b>	Lyophilized; PBS, no BSA, no azide, Purified: 1 mg/mL after reconstitution
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Storage Conditions</b>	Store at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Product origin</b>	Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
<b>Expected antibody applications</b>	IHC (frozen) (published for clone), Immunofluorescence (published for clone), Immuno-northern blot (published for clone), ELISA (published for clone), EMSA (published for clone), Flow (intracellular) (published for clone), IHC (FFPE) (published for clone)
<b>Shipping condition</b>	Room temperature