



Glowing products for science

# 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](#).

This datasheet was generated on September 1, 2025 at 12:10:16 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dsna-recombinant-monoclonal-mouse-antibody-rj2-lyophilized-biotium-choice/>

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

## Product attributes

Antibody number	#P025
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Clonality	Recombinant Monoclonal
Host species	Mouse
Clone	rJ2
Isotype	IgG2a, kappa
Antibody reactivity (target)	dsRNA
Synonyms	Double-stranded RNA; dsRNA
Antibody target cellular localization	Cytoplasm
Verified antibody applications	Dot blot (verified), IF (verified)
Positive control	poly(I-C) dot blot, HeLa cells
Antibody application notes	Stains endogenous dsRNA in HeLa cells by immunofluorescence. Selectively detects dsRNA on dot blot.
Antibody research areas	RNA Analysis & Detection, Vaccine Research, Virology
Antibody/conjugate formulation	Lyophilized; PBS, no BSA, no azide, Purified: 1 mg/mL after reconstitution
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Storage Conditions	Store at -35°C to -5°C
Regulatory status	For research use only (RUO)
Product origin	Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Expected antibody applications	IHC (frozen) (published for clone), Immunofluorescence (published for clone), ELISA (published for clone), EMSA (published for clone), Flow (intracellular) (published for clone), IHC (FFPE) (published for clone)
Filter by antibody application	Flow, intracellular (published), IHC, FFPE (published), IHC, frozen (published), Dot blot (verified), Immuno-northern blot (published), ELISA (published), EMSA (published), IF (verified)
Shipping condition	Room temperature