NeuN Recombinant Monoclonal Mouse Antibody (rA60) - Biotium Choice

NeuN (rA60) is a recombinant mouse monoclonal antibody that recognizes neuronal nuclear antigen protein NeuN. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

a fortun

Product Description

NeuN Recombinant Monoclonal Mouse Antibody (rA60) is a recombinant mouse monoclonal antibody that recognizes neuronal nuclear antigen protein NeuN. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to a selection of CF® Dyes.. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including near-IR CF®740
- Suitable for immunohistochemistry, immunofluorescence, and western

RNA binding protein fox-1 homolog 3 (RBFOX3), also known as NeuN, is a neuronal nuclear antigen encoded by the *Rbfox3* gene in mice. NeuN localizes to the nucleus and perinuclear cytoplasm of most CNS neurons, with notable absence in Cajal-Retzius, Purkinje, and photoreceptor cells. It is broadly expressed in brain regions (cortex, hippocampus, thalamus, striatum, cerebellum, spinal cord) from embryonic stages (E9.5–E12.5) through adulthood, and its expression is enhanced by retinoic acid. NeuN regulates alternative splicing, including *RBFOX2* isoforms targeted for nonsense-mediated decay. Initially identified by the monoclonal antibody A60, NeuN is now recognized as RBFOX3, and anti-NeuN antibodies label most neuronal populations. Knockout mice show reduced brain weight, impaired neurofilament expression, decreased white matter, seizure susceptibility, and reduced anxiety behaviors. Two isoforms (46 and 48 kDa) are present in nucleus and cytoplasm, with differing concentrations.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable CF® Dyes for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

Product attributes

Call us: 800-304-5357

Antibody number	P037
Antibody reactivity (target)	NeuN
Biotium Choice Antibody	Primary
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rA60
Isotype	lgG1
Molecular weight	50 kDa
Synonyms	Fox-1 homolog C, Fox-3, Hexaribonucleotide-binding protein 3, NeuN antigen, Neuronal nuclei antigen, RNA binding protein fox-1 homolog 3
Human gene symbol	RBFOX3
SwissProt	Q8BIF2
Immunogen	Purified nuclei from mouse brain
Antibody target cellular localization	Cytoplasm, Nucleus
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Ferret, Human, Mouse, Porcine, Rat, Salamander
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user
Positive control	Brain
Shipping condition	Room temperature
Storage Conditions	Store conjugates at 2 $^{\circ}\text{C}$ to 8 $^{\circ}\text{C},$ Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Antibody research areas	Neuroscience
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line
Cell/tissue expression	Brain, Neurons
Expected antibody applications	IHC (published for clone), IF (published for clone), WB (published for clone)

Email: btinfo@biotium.com

NeuN Recombinant Monoclonal Mouse Antibody (rA60) - Biotium Choice

Conjugation Ex/Em Conc. Size Catalog No. Dye Features

CF-05687 This datasheet was generated on November 2, 2025 at 10:17:14 AM. Visit product page to check for updated information before use.

Product link: https://biotium.com/product/neun-recombinant-monoclonal-mouse-antibody-ra60-biotium-choice/