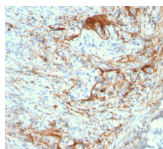


# Elastin Recombinant Monoclonal Rabbit Antibody (ELN/3131R)



## Product Description

Elastin is a polymeric protein and is the main component of the extracellular matrix of arteries. It is synthesized and secreted as a soluble, single-chain protein (tropoelastin), which undergoes a number of post-translational modifications prior to its organization into an elastic fiber in the extracellular space. Elastin performs a regulatory function during arterial development by controlling proliferation of smooth muscle and stabilizing arterial structure. It imparts elasticity to the connective tissue. During aging, the elasticity of connective tissue is reduced because of the cross-linking of collagenous fibers with elastin. The abnormal accumulation of elastic tissue in blood vessels is found in atherosclerosis and hypertension.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

**Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 3131, Anti-Elastin (ELN/3131R)**

## Product attributes

Antibody number	#3131
Antibody reactivity (target)	Elastin
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	ELN/3131R
Isotype	IgG
Molecular weight	70 kDa
Synonyms	Elastin; ELN; SVAS; Tropoelastin; WBS; WS
Human gene symbol	ELN
Entrez gene ID	2006
SwissProt	P15502
Unigene	647061
Immunogen	Recombinant full-length human Elastin protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Extracellular matrix
Species reactivity	Human
Positive control	A549 or U-2197 cells. Liver, Kidney, Lung or Intestine.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 13, 2025 at 12:47:43 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/elastin-recombinant-monoclonal-rabbit-antibody-eln-3131r/>