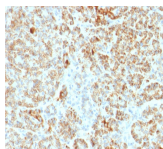


# Frataxin Monoclonal Mouse Antibody (FXN/2124)



## Product Description

Frataxin is a nuclear-encoded mitochondrial protein implicated in Friedreich's ataxia (FRDA), a human autosomal recessive neurodegenerative disease. Lack of Frataxin causes iron to accumulate in the mitochondrial matrix suggesting that Frataxin is involved in mitochondrial iron homeostasis and possibly in iron transport. Frataxin has an alpha-beta fold consisting of two helices flanking an antiparallel beta sheet.

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 2190, Anti-Faxin (FXN/2124)**

## Product attributes

Antibody number	#2190
Antibody reactivity (target)	Faxin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FXN/2124
Isotype	IgG2b, kappa
Molecular weight	14 kDa
Synonyms	CyaY; d-FXN; FARR; Frataxin mature form; Frataxin (81-210); FRDA; Friedreich ataxia protein; Fxn; l-FXN; m56-FXN; m78-FXN; m81-FXN; X25
Human gene symbol	FXN
Entrez gene ID	2395
SwissProt	Q16595
Unigene	20685
Immunogen	Recombinant fragment of human Frataxin (FXN) protein (around aa 57-210) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Mitochondria
Species reactivity	Human
Positive control	HepG2, Panc-2, HL-60 or HEK293 cells. Pancreas and Liver.
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
Validated in protein array	Monospecific
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 8, 2025 at 01:43:36 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/frataxin-monoclonal-mouse-antibody-fxn-2124/>