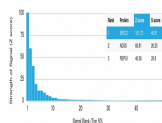




Glowing products for science

ERCC1 / RAD10 Monoclonal Mouse Antibody (ERCC1/2318)



Product Description

Recognizes a protein of 110 kDa, identified as Excision Repair Cross Complementing 1 (ERCC1). It is a mammalian nucleotide excision repair (NER) enzyme involved in repair of damaged DNA. ERCC1 is a homologous to RAD10 in *Saccharomyces cerevisiae*, which is required in mitotic intrachromosomal recombination and repair. ERCC1 is required in repair of cisplatin-induced DNA adducts and ultraviolet (UV)-induced DNA damage. High expression of ERCC1 has been linked to tumor progression in a variety of cancers including non-small cell lung cancer (NSCLC), squamous cell carcinoma of the head, ovarian cancer and esophageal cancer.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2318, Anti-ERCC1|RAD10 (ERCC1/2318)

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

Product attributes

Antibody number	#2318
Antibody reactivity (target)	ERCC1, RAD10
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ERCC1/2318
Isotype	IgG2b, kappa
Molecular weight	~110 kDa
Synonyms	COFS4; DNA excision repair protein ERCC1; ERCC1; Excision repair cross complementing 1; RAD10; UV20
Human gene symbol	ERCC1
Entrez gene ID	2067
SwissProt	P07992
Unigene	435981
Immunogen	Recombinant fragment (around aa 191-281) of human ERCC1 protein (exact sequence is proprietary)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	HeLa, HepG2 cells. Tonsil or Prostate.
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 (<i>Bos taurus</i>), 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	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
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 January 8, 2026 at 03:23:22 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ercc1-rad10-monoclonal-mouse-antibody-ercc1-2318/>