## Cathepsin D Monoclonal Mouse Antibody (CTSD/3083)



## **Product Description**

Cathepsin D is a ubiquitously expressed lysosomal aspartyl protease involved in the normal degradation of proteins. It is synthesized as an inactive 43 kDa preprocathepsin D that is cleaved and glycosylated to form a 46 kDa procathepsin D and then further cleaved to produce 28 kDa and 15 kDa subunits (heavy and light chains, respectively). Cathepsin D exhibits pepsin-like activity and plays a role in protein turnover and in the proteolytic activation of hormones and growth factors. Mutations in this gene play a causal role in neuronal ceroid lipofuscinosis-10 and may be involved in the pathogenesis of several other diseases, including breast cancer and possibly Alzheimer's disease.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3083, Anti-Cathepsin D (CTSD/3083)

## Product attributes

Call us: 800-304-5357

Froduct attributes		
Antibody number	#3083	
Antibody reactivity (target)	Cathepsin D	
Antibody type	Primary	
Host species	Mouse	
Clonality	Monoclonal	
Clone	CTSD/3083	
Isotype	IgG2b, kappa	
Molecular weight	46 kDa (Procathepsin D); 28 kDa (Cathepsin D)	
Synonyms	Cathepsin D; CTSD; Lysosomal aspartyl peptidase	
Human gene symbol	CTSD	
Entrez gene ID	1509	
SwissProt	P07339	
Unigene	654447	
Immunogen	Recombinant fragment of human Cathepsin D protein (around as 104-250) (exact sequence is proprietary)	
Verified antibody applications	IHC (FFPE) (verified), WB (verified)	
Antibody target cellular localization	Secreted (extracellular)	
Species reactivity	Human	
Positive control	K-562, A431 or SK-BR3 cells. Kidney, Liver, Lung or Breast.	
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.	

Email: btinfo@biotium.com

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 November 2, 2025 at 03:51:39 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cathepsin-d-monoclonal-mouse-antibody-ctsd-3083/">https://biotium.com/product/cathepsin-d-monoclonal-mouse-antibody-ctsd-3083/</a>