

Luteinizing Hormone / Choriogonadotropin Receptor Monoclonal Mouse Antibody (LHCGR/1416)

Product Description

This antibody recognizes a protein of 78-85 kDa, which is identified as luteinizing hormone/choriogonadotropin receptor. Luteinizing hormone plays a role in spermatogenesis and ovulation by stimulating the testes and ovaries to produce steroids. Choriogonadotropin production in the placenta maintains estrogen and progesterone levels during the first trimester of pregnancy. Ovaries and testes abundantly express luteinizing hormone/choriogonadotropin receptor (LHCGR) as a seven transmembrane, G protein-coupled receptor glycoprotein. LHCGR influences the protective effect of pregnancy and Gonadotropin against breast cancer. The expression of LHCGR on breast carcinoma correlates in part to the degree of tumor differentiation. LHCGR -positive breast tumors occur more frequently in tumors with greater cell differentiation in premenopausal women.

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 1416, Anti-Choriogonadotropin Receptor (LHCGR/1416)

Product attributes

Antibody number	#1416
Antibody reactivity (target)	Choriogonadotropin Receptor, Luteinizing Hormone
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LHCGR/1416
Isotype	IgG1, kappa
Molecular weight of antigen	78-85 kDa
Synonyms	Gonadotropin receptor; GTHR-II; HHG; LCGR; LGR2; LH-R; LH/CG-R; LH/CGR; LHCGR; LHR; LHRHR; LSH-R; Luteinizing hormone receptor; Luteinizing hormone/choriogonadotropin receptor; Lutropin choriogonadotropic receptor; Lutropin-choriogonadotropic hormone receptor; ULG5
Human gene symbol	LHCGR
Entrez gene ID	3973
SwissProt	P22888
Unigene	468490
Immunogen	Purified recombinant N-terminal fragment of human LHCGR (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane
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	Ovary or testis
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
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: 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
Antibody research areas	Cancer, Endocrinology
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.
Tumor expression	Breast cancer

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 June 19, 2026 at 01:01:38 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mouse-monoclonal-anti-luteinizing-hormone-choriogonadotropin-receptor-lhgr1416/>