Luteinizing Hormone / Choriogonadotropin Receptor Monoclonal Mouse Antibody (LHCGR/1417)

Product Description

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 1417, Anti-Choriogonadotropin Receptor (LHCGR/1417)

Product attributes

Tumor expression

| Product attributes | | | | |
|---------------------------------------|---|--|--|--|
| Antibody number | #1417 | | | |
| Antibody reactivity (target) | Choriogonadotropin Receptor, Luteinizing Hormone | | | |
| Antibody type | Primary | | | |
| Host species | Mouse | | | |
| Clonality | Monoclonal | | | |
| Clone | LHCGR/1417 | | | |
| Isotype | IgG1, kappa | | | |
| Molecular weight | 78-85 kDa | | | |
| Synonyms | Gonadotropin receptor; GTHR-II; HHG; LCGR; LGR2; LH-R; LH/GGR; LH/CGR; 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 | 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 i 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. | | | |

Breast cancer

Call us: 800-304-5357 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 5, 2025 at 11:42:37 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mouse-monoclonal-anti-luteinizing-hormone-choriogonadotropin-receptor-lhcgr1417/