

# FSH Receptor Monoclonal Mouse Antibody (FSHR/1400)

## Product Description

Follicle-stimulating hormone receptor (FSHR) is a 695 amino acid G protein coupled receptor. FSH binds to the receptor in a hand-clasp fashion via its alpha and beta subunits. While the beta subunit of FSH is involved in the binding of FSH to the receptor, the alpha subunit stabilizes this interaction. Linkage studies suggest that a missense mutation in the FSHR gene can cause reduced FSH binding affinity and lead to a condition known as hypergonadotropic ovarian dysgenesis (ODG). In males however, this mutation does not appear to have a detrimental affect on fertility. It is believed that a mutation in the FSHR gene is also associated with ovarian hyperstimulation syndrome; a condition characterized by the presence of multiple serous and hemorrhagic follicular cysts lined by luteinized cells. 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 1400, Anti-FSH Receptor (FSHR/1400)**

## Product attributes

Antibody number	#1400
Antibody reactivity (target)	FSH Receptor
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FSHR/1400
Isotype	IgG1
Molecular weight	75 kDa
Synonyms	Follicle-stimulating hormone receptor; Folitropin receptor; FSH receptor; FSH-R; FSHRO; LGR1; ODG1; ovarian dysgenesis 1
Human gene symbol	FSHR
Entrez gene ID	2492
SwissProt	P23945
Unigene	1428
Immunogen	Recombinant human full-length FSHR protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology (formalin) 1-2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	Uterine Carcinoma
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	Endocrinology

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>
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, LI-COR Bioscience.