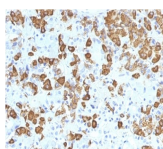


# Growth Hormone Monoclonal Mouse Antibody (GH/1371)



## Product Description

Pituitary growth hormone (GH) plays a crucial role in stimulating and controlling the growth, metabolism and differentiation of many mammalian cell types by modulating the synthesis of multiple mRNA species. These effects are mediated by the binding of GH to its membrane-bound receptor, GHR, and involve a phosphorylation cascade that results in the modulation of numerous signaling pathways. GH is synthesized by acidophilic or somatotrophic cells of the anterior pituitary gland. Anti-GH is a useful marker in classification of pituitary tumors and the study of pituitary disease (acromegaly).

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 1371, Anti-Growth Hormone (GH/1371)**

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

## Product attributes

|  |  |
|--|--|
| <b>Antibody number</b>                       | #1371  |
| <b>Antibody reactivity (target)</b>          | Growth Hormone   |
| <b>Antibody type</b>                         | Primary  |
| <b>Host species</b>                          | Mouse  |
| <b>Clonality</b>                             | Monoclonal   |
| <b>Clone</b>                                 | GH/1371  |
| <b>Isotype</b>                               | IgG2b, kappa   |
| <b>Molecular weight of antigen</b>           | 20 kDa   |
| <b>Synonyms</b>                              | GH; GH-N; GH1; GHN; Growth hormone 1; Growth hormone; Growth hormone; pituitary; HG1; hGH-N; IGHD1B; Pituitary growth hormone; RNGTHGP; Somatotropin   |
| <b>Human gene symbol</b>                     | GH1  |
| <b>Entrez gene ID</b>                        | 2688   |
| <b>SwissProt</b>                             | P01241   |
| <b>Unigene</b>                               | 655229   |
| <b>Immunogen</b>                             | A recombinant human Growth Hormone (GH) fragment (aa58-187) (exact sequence is proprietary)  |
| <b>Antibody target cellular localization</b> | Secreted (extracellular)   |
| <b>Verified antibody applications</b>        | IHC (FFPE) (verified)  |
| <b>Species reactivity</b>                    | Human  |
| <b>Antibody application notes</b>            | 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): 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, 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 |
| <b>Positive control</b>                      | Pituitary cells or Pituitary   |
| <b>Shipping condition</b>                    | Room temperature   |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C  |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended  |
| <b>Regulatory status</b>                     | For research use only (RUO)  |
| <b>Antibody/conjugate formulation</b>        | 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  |
| <b>Validated in protein array</b>            | Monospecific   |
| <b>Antibody research areas</b>               | Endocrinology  |
| <b>Cell/tissue expression</b>                | Pituitary gland  |
| <b>Product origin</b>                        | 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 | <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>  |
| BNC74             | CF®740             | 742/767    | 633-685    | 775/50                   | <a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on June 14, 2026 at 12:13:55 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mouse-monoclonal-anti-growth-hormone-gh1371/>