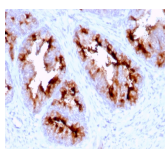


FOLH1 / PSMA Monoclonal Mouse Antibody (FOLH1/2363)



Product Description

Folate hydrolase 1 (FOLH1), also known as Prostate-specific membrane antigen (PSMA), is a type II transmembrane glycoprotein belonging to the M28 peptidase family. FOLH1 has two enzymatic activities, one as a prostate-specific integral membrane folate hydrolase and the other as a carboxypeptidase. In the prostate the protein is up-regulated in cancerous cells and is used as an effective diagnostic and prognostic indicator of prostate cancer.

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 2363, Anti-FOLH1|PSMA (FOLH1/2363)

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

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #2363 |
| Antibody reactivity (target) | FOLH1, PSMA |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | FOLH1/2363 |
| Isotype | IgG2b, kappa |
| Molecular weight of antigen | 100 kDa |
| Synonyms | Cell growth-inhibiting gene 27 protein (GIG27); Folate hydrolase 1 (FOLH1); Foly/poly-gamma-glutamate carboxypeptidase (FGCP); Glutamate carboxylase II (GCP1I); Glutamate carboxypeptidase 2 (GCP2); Glutamate carboxypeptidase II Membrane glutamate carboxypeptidase N-acetylated-alpha-linked acidic dipeptidase I (NAALAD1 or NAALADase); Prostate-specific membrane antigen (PSM or PSMA); Pteroyl/poly-gamma-glutamate carboxypeptidase |
| Human gene symbol | FOLH1 |
| Entrez gene ID | 2346 |
| SwissProt | Q04609 |
| Unigene | 654487 |
| Immunogen | Recombinant human FOLH1 protein fragment (around aa 232-433) (exact sequence is proprietary) |
| Verified antibody applications | IHC (FFPE) (verified), WB (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | LNCap or HepG2 cells. Prostate 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 |
| 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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Cell/tissue expression | Prostate |
| Tumor expression | Prostate 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 08:32:54 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/folh1-psma-monoclonal-mouse-antibody-folh1-2363/>