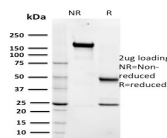


## Maltose Binding Protein / MBP-probe Monoclonal Mouse Antibody (R29.6)



### Product Description

Plasmid vectors for the expression of coding regions of eukaryotic genes in bacterial, insect and mammalian hosts are in common usage; such expression vectors frequently encode hybrid fusion proteins consisting in part of prokaryotic and in part, eukaryotic specified proteins. One such system utilizes maltose binding protein (MBP), the 370 amino acid product of the E. coli mal E gene. Plasmid vectors have been constructed utilizing the MBP domain that allow the synthesis of high levels of MBP-fusion proteins that can be Purified in a one step procedure by affinity chromatography crosslinked amylose resin. Once bound to amylose, the MBP protein can then be separated from the target protein by cleavage by coagulation factor Xa at a specific four residue site. Alternatively, the intact fusion protein can be specifically eluted from the resin by the addition of excess free maltose. Subsequent to elution, MBP fusion protein can be visualized either by western blot analysis or immunoprecipitation using antibodies specific for the MBP-tag. Expression systems utilizing the MBP fusion tag include pCG-806fx and pMal vectors. 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.

### Product attributes

|                                       |   |
|---------------------------------------|---|
| Antibody number                       | #2847   |
| Antibody reactivity (target)          | Maltose Binding Protein, MBP-probe  |
| Antibody type                         | Primary   |
| Host species                          | Mouse   |
| Clonality                             | Monoclonal  |
| Clone                                 | R29.6   |
| Isotype                               | IgG1, kappa   |
| Molecular weight                      | 40 kDa  |
| Synonyms                              | ECK4026; JW3994; MalE; malJ; Maltodextrin binding protein; Maltose ABC transporter periplasmic protein; Maltose binding periplasmic protein; Periplasmic maltose binding protein  |
| Human gene symbol                     | Not Known   |
| Entrez gene ID                        | Not Known   |
| SwissProt                             | Not Known   |
| Unigene                               | Not Known   |
| Immunogen                             | MOS maltose binding protein fusion protein.   |
| Verified antibody applications        | WB (verified)   |
| Antibody target cellular localization | Periplasm   |
| Antibody application notes            | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL; Western blot 0.5-1.0 ug/mL; Optimal dilution for a specific application should be determined. |
| Positive control                      | Saas-2 cells. MBP fusion proteins.  |
| Shipping condition                    | Room temperature  |
| Storage Conditions                    | Note: store BSA-free antibodies at -10 to -35°C; Store at 2 to 8 °C; Protect fluorescent conjugates from light  |
| 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   |
| Shelf life                            | Guaranteed for at least 24 months from date of receipt when stored as recommended   |

| 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.

This datasheet was generated on March 29, 2024 at 03:24:33 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/maltose-binding-protein-mbp-probe-monoclonal-mouse-antibody-r29-6/>