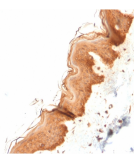


Cytokeratin 4 Monoclonal Mouse Antibody (KRT4/2804)



Product Description

Cytokeratin 4 (KRT4) is an intermediate filament protein associated with Cytokeratin 13 (KRT13). It is expressed in suprabasal cells of non-keratinized stratified squamous epithelium of esophagus, cornea, anus, larynx, pharynx and tongue. Decreased expression of KRT4 is associated with head and neck squamous carcinoma. It is helpful in differentiation of squamous cell carcinoma of esophagus origin from that of thyroid origin.

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 2804, Anti-Cytokeratin 4 (KRT4/2804)

Product attributes

| | |
|--|---|
| Antibody number | #2804 |
| Antibody reactivity (target) | Cytokeratin 4 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | KRT4/2804 |
| Isotype | IgG1, kappa |
| Molecular weight of antigen | 67 kDa |
| Synonyms | CK4; CYK4; Cytokeratin 4; Keratin 4; Keratin type II cytoskeletal 4; KRT4 |
| Human gene symbol | KRT4 |
| Entrez gene ID | 3851 |
| SwissProt | P19013 |
| Unigene | 654610; 706527 |
| Immunogen | Recombinant fragment (around aa 181-292) of human KRT4 protein (exact sequence is proprietary) |
| Verified antibody applications | Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | HeLa, A549 or A431 cells. Skin, Cervix or Esophagus. |
| 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 | Epithelial cells, Eye |
| Antibody research areas | Cancer, Cytoskeleton |
| Tumor expression | Squamous cell carcinoma |

| 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 5, 2026 at 01:22:09 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cytokeratin-4-monoclonal-mouse-antibody-krt4-2804/>