

XPB ((Xeroderma pigmentosum group B) Monoclonal Antibody

Cat No: G01069 - 100 μ L

General Data

| | |
|---------------------|---|
| Shipping: | wet ice |
| Stability: | Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative |
| Formulation: | Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles |
| Host: | Mouse |
| Antigen: | Recombinant human TFIIHp89/XPB protein. |
| Clone: | 1B3 |
| Isotype: | IgG1, κ |



| | |
|------------------------|-------------------------------------|
| Application(s): | ELISA |
| | Western Blot |
| | Immunoprecipitation |
| | Immunocytochemistry |
| | Immunohistochemistry |
| | ChIP |
| | Recommended dilutions: 1/500-1/5000 |

Specificity: human TFIIHp89/XPB.

Scientific Literature

Schaeffer L, Roy R, Humbert S, Moncollin V, Vermeulen W, Hoeijmakers JH, Chambon P, Egly JM. DNA repair helicase: a component of BTF2 (TFIIH) basic transcription factor. Science. 1993 Apr 2;260(5104):58-63

Le May N, Mota-Fernandes D, Vélez-Cruz R, Iltis I

FP/08/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>