

BRM (human), (recombinant)

BRM is a subunit of SWI-SNF complex, an essential effector of chromatin architecture and transcriptional regulation. It belongs to a family of SNF2-related nucleosome remodeling enzymes (DNA-dependent ATPases). These ATP-dependent motor proteins are major factors of chromatin structure and transcriptional control in vivo. SNF2-related factors have been implicated in a variety of cancers in humans.

Ordering Information

[Order Online »](#)

| | |
|-----------------|------|
| ENZ-PRT334-0050 | 50µl |
|-----------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------------|---|
| Handling | Thaw and freeze quickly, in room-temperature water bath and liquid nitrogen, respectively; keep on ice when not in storage; split in small aliquots upon first thawing; avoid excessive freeze-thaw cycles. |
| Long Term Storage | -80°C |
| Shipping | Dry Ice |

Product Details

| | |
|-------------------------|---|
| Alternative Name | Probable global transcription activator SNF2L2, BAF190B, SMCA2, SNF2- α , SNF2A, SNF2LA, SMARCA2 |
| Formulation | Liquid. In 20 mM Tris-HCl, pH 7.9, 150 mM NaCl, 15% glycerol, 2 mM MgCl ₂ , 0.2 mM EDTA, 1 mM DTT, 10 mM beta-glycerophosphate, 0.4 mM PMSF, 1 mM benzamidine and 0.4 mg/ml recombinant human insulin. |
| MW | ~181 kDa |
| Purity Detail | Purified by immunoaffinity chromatography. |
| Source | Produced in insect (Sf9) cells. Human BRM (aa 1-1572) fused to 6xHis and FLAG-tag at the N-terminus. |
| UniProt ID | P51531-2 |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com