

BLM Helicase, Active, Human

Part number: 2353 (5 µg), 2354 (50 µg)
 Lot number: 14975

BLM Helicase, Active, Human

General: BLM Helicase is a human enzyme with a C-terminal 6xHis tag.

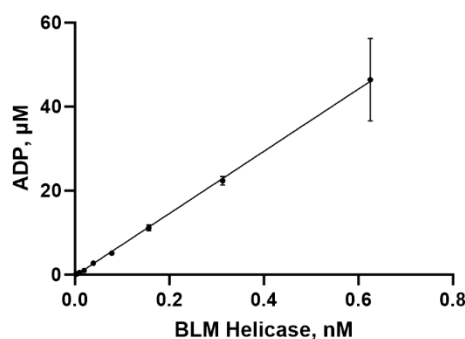
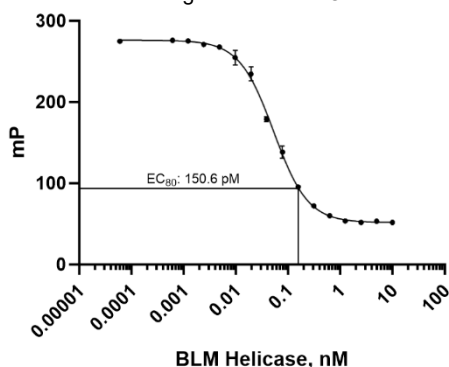
Storage and Handling: Store at -80°C. Product should remain active for 6 months after purchase date with minimal freeze-thaws.

Purity: ≥85% as determined by SDS-PAGE

Expression: Sf9 (baculovirus)

Description	Composition	Concentration	Rate*	Storage
BLM Helicase, 76.16 kDa	50 mM Tris-HCl, 500 mM NaCl, 5% Glycerol, pH 8.0	0.3 mg/mL (3.94 µM)	20.2 sec ⁻¹	-80°C

*Rate of ADP formation measured using Transcreener® ADP FP Assay



BLM Helicase, Human: Functional Activity

Activity was measured using the Transcreener® ADP² FP Assay (100 µM ATP and 1 µg/mL Sheared Salmon Sperm DNA in 1X Enzyme Assay Buffer H (50 mM Tris, 5 mM MgCl₂, 100 mM NaCl and 0.01% Triton X-100, pH 8.0) with 1 mM DTT; 10 µL Enzyme Reaction incubated @ 30°C for 60 minutes in a Corning® low-volume 384-well plate (#4514). 20 µL Complete Reaction read on CLARIOstar Plus instrument. Fluorescent polarization values were converted to ADP formation with a 100 µM ADP Standard Curve and ADP formation was plotted using Prism (GraphPad). Under these conditions, it is recommended to use BLM Helicase in the 10 µL Enzyme Reaction at a concentration between the EC₅₀ and EC₈₀ of the enzyme titration for a robust signal in the linear range.

BLM Helicase, Human: Amino Acid Sequence

Number of Amino Acids: 668 (aa 630-1290)

MSAQNLASRNKHERFQSLSPHTKEMMKIFHKKFGLHNFRTNQLEAINAALLGEDCFILMPTGGGKSLCYQLPACVSPGVTVVISPLRSLIVD
 QVQKLTSLDIPATYLTGDKTDSEATNIYLQLSKKDPIIKLLYTPEKICASNRLISTLENLYERKLLARFVIDEAHCVSQWGHDFRQDYKRMNML
 RQKFPSPVPMALTATANPRVQKDILTQLKILRPQVFSMSFNHNLKYYVLPKKPKKVAFDCLWIRKHHYPYDSGIIYCLSRRECDTMDATLQR
 DGLAALAYHAGLSDSARDEVQKQWVNDGCGVICATIAFGMGIDKPDVRFVIHASLPKSVGGYQESGRAGRDEISHCLLFYTYHDVTRLKR
 LIMMEKDGNNHHTRETFNNLYSMVHYCENTECRRILLAYFGENGFNPDFCKKHPDVSCDNCCKTKDYKTRDVTDDVKISIVRFVQEHSSSQ
 GMRNIKHVGPGRFTMNLVDIFLGSKSAKIQSGIFGKGSAYSRRHAERLFKKLILDKILDEDLYINANDQAIAYVMLGNKAQTVLNGNLKVDF
 METENSSSVKKQKALVAKVSQREEMVKKCLGELTEVCKSLGKVGHVHYFNIFNTVTLKLAESLSDPEVLLQIDGVTEDEKLEKYGAEIVSVL
 QKYSEWTSHHHHH

User Notification

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. BellBrook Labs LLC will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use, or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, BellBrook Labs LLC is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, BellBrook Labs LLC, 1232 Fourier Drive, Suite 115, Madison, Wisconsin 53717. Phone (608)443-2400. Fax (608)441-2967.

©2025 BellBrook Labs. All rights reserved.