

Rbx1 (human), (recombinant) (His- tag)

Cullin-based E3 ligases, such as SCF (Skp1-Cdc53-F-box protein), are a large family of multi-subunit ubiquitin ligases with diverse cellular functions, including the regulation of the cell cycle, of the DNA damage response, and of various transcription factors. These ligases are composed of one of six mammalian cullin homologs (Cul1, Cul2, Cul3, Cul4a, Cul4b, and Cul5), the Ring finger containing protein Rbx1/Roc1, and cullin homolog-specific adaptor and substrate recognition subunits. Activation of cullin-based ligases require the covalent modification of a conserved lysine residue in the cullin protein with the ubiquitin-like protein NEDD8. This modification is catalyzed by the Rbx1 subunit.

Ordering Information

[Order Online »](#)

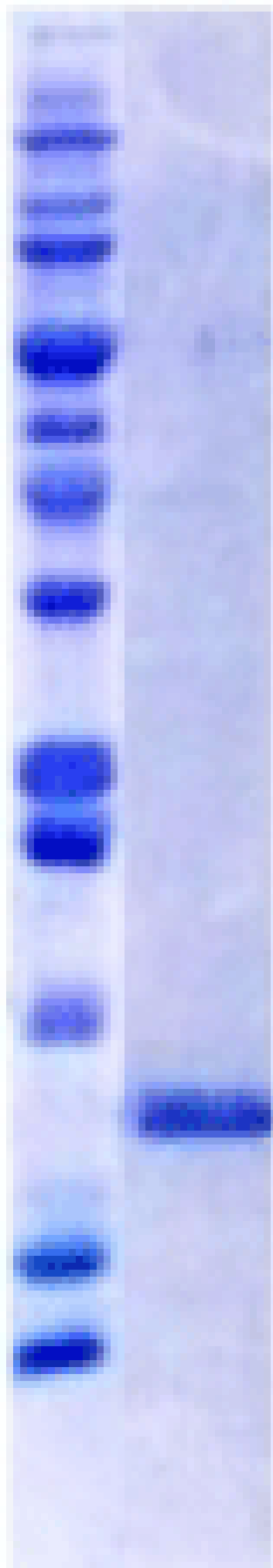
BML-UW0395-0025	25µg
------------------------	------

Manuals, SDS & CofA

[View Online »](#)

1

2



Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ring box protein-1, ROC1, Regulator of cullins 1, Ring finger protein 75
Formulation	Liquid. In 50mM TRIS, pH 8.0, containing 50mM NaCl and 1mM DTT.
MW	~14kDa
Purity	≥98% (SDS-PAGE)
Purity Detail	Purified by nickel sepharose affinity chromatography.
Source	Produced in <i>E. coli</i> BL21 (λDE3).
UniProt ID	P62877

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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