

UbcH10 (human), (recombinant) (untagged)

The anaphase-promoting complex (APC) coordinates mitosis by sequentially promoting the degradation of key cell-cycle regulators *via* the ubiquitin-proteasome pathway, utilizing the ubiquitin E2 conjugating enzyme UbcH10 and K11-linked polyubiquitin chain modification of substrate proteins. Following the degradation of its substrates, the APC catalyzes the autoubiquitination of UbcH10 resulting in its degradation and subsequent inactivation of the APC.

Aberrant performance or modulated cellular levels of UbcH10 or any of its associated components is likely to have significant consequences for cell cycle regulation.

UbcH10 expression has been shown to be a valid indicator of the proliferative and aggressive status of thyroid carcinomas and its overexpression is thought to play a role ovarian carcinogenesis. In contrast expression levels of UbcH10 are extremely low in many normal cells.

Such observations suggest the possibility of using UbcH10 detection as a marker for the diagnosis and prognosis of neoplastic diseases and open the possibility of a therapy of some carcinomas based on the suppression of the UbcH10 synthesis and/or function.

Citations: 2

[View Online »](#)

Ordering Information

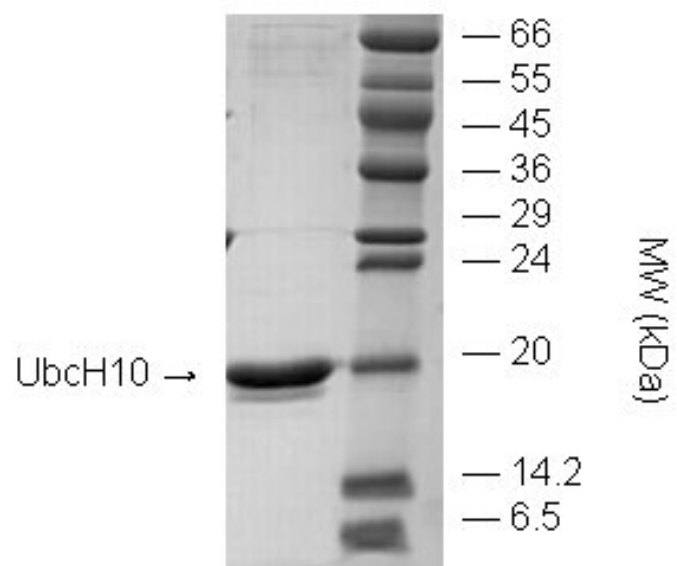
[Order Online »](#)

BML-UW0960-0100

100µg

Manuals, SDS & CofA

[View Online »](#)



Coomassie stained gel of UbchH10 (5µg, BML-UW0960)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored, as supplied, 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	UBE2C, Ubiquitin-protein ligase C
Application Notes	E2 conjugating enzyme for ubiquitinylation reactions.
Formulation	Liquid. In PBS.
MW	~19.7kDa
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	O00762



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)