

UBE2K, His₆

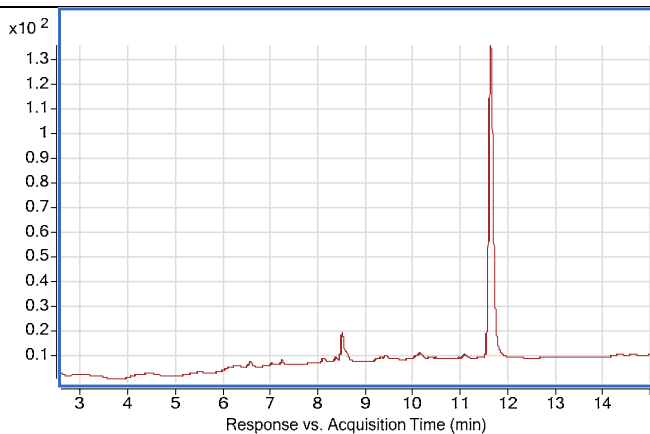
Cat. # UB204H

Background: UBE2K, also known as UbcH1 or E2-25K, is an E2 (ubiquitin conjugating enzyme) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E3 enzymes. Once an E1 is carrying an activated ubiquitin in a ubiquitin-E1 thiolester complex, this activated ubiquitin is transferred to the active site cysteine of the E2 to form a ubiquitin-E2 thiolester complex. With the aid of an E3, the activated ubiquitin is then targeted to a specific substrate lysine. UBE2K is a biochemically important E2 due to its ability to form free K48-linked polyubiquitin chains in solution in the absence of an E3. This ability is attributed to UbcH1 possessing a C-terminal UBA domain, in addition to its E2 catalytic domain.

Application: Ubiquitin ligation reactions

Product Information

Source:	Human, recombinant; Accession No. P61086
Organism:	<i>E. coli</i>
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	23,400.7 Da
Tag	His ₆
Physical State:	Liquid,
Quantity:	20 µL or 75 µL of a 40 µM solution (0.8 and 3 nmoles, respectively)
Solubility:	>3 mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles



All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved