

UBE2K

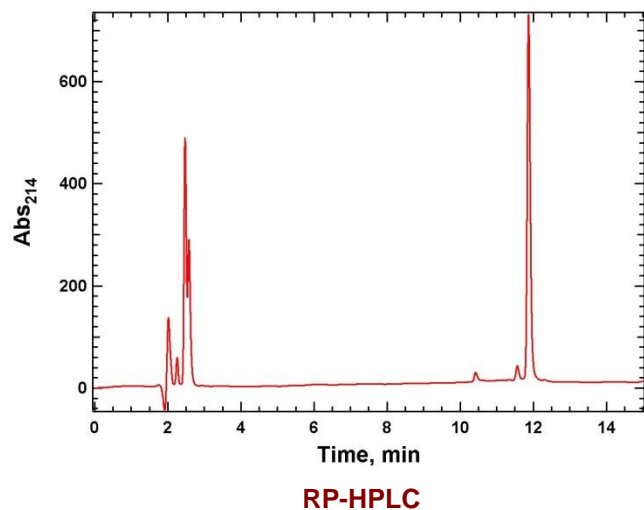
Cat. # UB204

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

Organism	Human, recombinant; Accession No. P61086
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	22,406.7 Da
Tag	none
Physical State:	Liquid, TBS
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