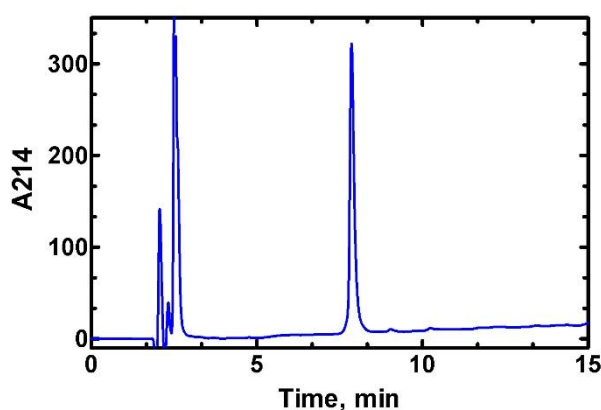


Background	UBE2L6 (also known as Ubch8) is an E2 ubiquitin-conjugating enzyme involved in the conjugation of ubiquitin to target substrates, in collaboration with E1 and E3 enzymes. UBE2L6 is also responsible for the conjugation of the ubiquitin-like protein (Ubl) ISG15 to its substrates (a process known as ISGylation), in concert with the ISG15 E1 activating enzyme. ISG15 is an interferon-induced Ubl that plays a role in innate immunity pathways.
Alternate Names	Ubch8, Ubiquitin/ISG15-Conjugating Enzyme E2 L6, E2 Ubiquitin-Conjugating Enzyme L6, RIG-B, Retinoic Acid Induced Gene B Protein
Application(s)	Ubiquitin ligation reactions

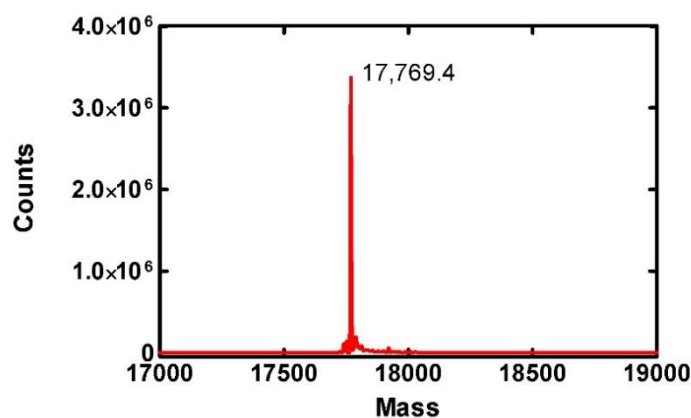
Product Specifications

Tag	None
Purity	≥ 90% by RP-HPLC
Molecular Weight	17,768.8 Da
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	Human, Accession No. O14933
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
Solubility	> 3 mg/ml
Stability & Storage	1 year at -80°C. Avoid repeated freeze/thaw cycles

Product QC



RP-HPLC



Deconvoluted mass spectrum

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved

References

1. Shi, H. X., Yang, K., Liu, X., Liu, X. Y., Wei, B., Shan, Y. F., Zhu, L. H., & Wang, C. (2024). HERC5-catalyzed ISGylation potentiates cGAS-mediated innate immunity. *Cell Reports*, 39(1), 110658.
2. Durfee, L. A., Kelley, M. L., & Huibregtse, J. M. (2008). The basis for selective E1–E2 interactions in the ISG15 conjugation system. *Journal of Biological Chemistry*, 283(35), 23895–23902.
3. Dastur, A., Beaudenon, S., Kelley, M., Krug, R. M., & Huibregtse, J. M. (2006). Herc5, an interferon-induced HECT E3 enzyme, is required for conjugation of ISG15 in human cells. *Journal of Biological Chemistry*, 281(7), 4334–4338.
4. Wong, J. J., Pung, Y. F., Sze, N. S., & Chin, K. C. (2006). HERC5 is an IFN-induced HECT-type E3 protein ligase that mediates type I IFN-induced ISGylation of protein targets. *Proceedings of the National Academy of Sciences*, 103(28), 10735–10740

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved