

Otulin (FAM105B)

Cat. # DB206

Background: Otulin was recently described as a deubiquitylase exhibiting high specificity for linear ubiquitin chains (M1)^(1,2). Specific interactions between the proximal ubiquitin and Otulin leads to activation of the catalytic triad. Otulin binds to the HOIP (RNF31) component of the linear chain synthesizing E3 ligase complex, LUBAC, and regulates its activity thus affecting NFκB regulated pathways and proinflammatory responses^(1,3). LUBAC regulation of NFκB is also involved in angiogenesis⁽²⁾.

Application: Specific cleavage of linear (M1) ubiquitin chains

Product Information

| | |
|--------------------------|---|
| Origin: | Human, recombinant. Accession number NP_612357 |
| Other names: | Gumby, FAM105B |
| Purity: | ≥ 95% by |
| Molecular Weight: | 40,579 Da |
| Physical State: | Liquid, 20mM Hepes, pH 7.5, 100mM NaCl, 5mM 2-mercaptoethanol |
| Quantity: | 25 µg |
| Solubility: | >20mg/mL |
| Storage: | -80° C. Avoid repeated freeze/thaw cycles |

References

1. Keusekotten, K., Elliott, P. R., Glockner, L., Fiil, B. K., Damgaard, R. B., Kulathu, Y., Wauer, T., Hospenthal, M. K., Gyrd-Hansen, M., Krappmann, D., Hofmann, K., and Komander, D. (2013) *Cell* **153**, 1312-1326
2. Rivkin, E., Almeida, S. M., Ceccarelli, D. F., Juang, Y. C., MacLean, T. A., Srikumar, T., Huang, H., Dunham, W. H., Fukumura, R., Xie, G., Gondo, Y., Raught, B., Gingras, A. C., Sicheri, F., and Cordes, S. P. (2013) *Nature* **498**, 318-324
3. Fiil, B. K., Damgaard, R. B., Wagner, S. A., Keusekotten, K., Fritsch, M., Bekker-Jensen, S., Mailand, N., Choudhary, C., Komander, D., and Gyrd-Hansen, M. (2013) *Mol Cell* **50**, 818-830

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2014 LifeSensors, Inc. All Rights Reserved