

## PLP2 (Papain like protease 2)

Cat. # DB603

**Background:** To date, five human coronaviruses (CoV) have been identified including the common human respiratory pathogen NL63, that has been shown to be associated with both upper and lower respiratory tract disease such as bronchiolitis and croup<sup>1-3</sup>. The prevalence of NL63 infections is illustrated by epidemiological studies reporting that NL63 CoV was detected in 0.3-10% of clinical specimens from both adults and infants diagnosed with respiratory tract diseases<sup>4-12</sup>. The NL63 genome encodes two papain like proteases, PLP1 and PLP2. PLP2 is a dual specificity isopeptidase with both deubiquitylase and deISGylase activity<sup>13,14</sup>. PLP2 cleaves Ub-PLA<sub>2</sub> and ISG15-PLA<sub>2</sub> with high and moderate efficiency respectively.

**Alternate names:** A4, A4-LSB, Differentiation-dependent protein A4, Intestinal membrane A4 protein, MGC126187, Proteolipid protein 2

### Product Information

<b>Molecular Weight:</b>	36.7kDa
<b>Quantity:</b>	25µg
<b>Physical State:</b>	Liquid
<b>Source:</b>	NL63-CoV Recombinant
<b>Tag:</b>	His6
<b>Activity:</b>	This enzyme is active in the Ub-CHOP and ISG15-CHOP assay.
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

### References

- 1) Van der Hoek, L., et al., *Identification of a new human coronavirus*. Nat Med, 2004. 10(4): p. 368-73.
- 2) Van der Hoek, L., et al., *Croup is associated with the novel coronavirus NL63*. PLoS Med, 2005. 2(8): p. e240.
- 3) Fouchier, R.A., et al., *A previously undescribed coronavirus associated with respiratory disease in humans*. Proc Natl Acad Sci U S A, 2004. 101(16): p. 6212-6.
- 4) Vabret, A., et al., *Human coronavirus NL63, France*. Emerg Infect Dis, 2005. 11(8): p. 1225-9.
- 5) Suzuki, A., et al., *Detection of human coronavirus-NL63 in children in Japan*. Pediatr Infect Dis J, 2005. 24(7): p. 645-6.
- 6) Kaiser, L., et al., *Human coronavirus NL63 associated with lower respiratory tract symptoms in early life*. Pediatr Infect Dis J, 2005. 24(11): p. 1015-7.
- 7) Ebihara, T., et al., *Detection of human coronavirus NL63 in young children with bronchiolitis*. J Med Virol, 2005. 75(3): p. 463-5.
- 8) Choi, E.H., et al., *The association of newly identified respiratory viruses with lower respiratory tract infections in Korean children, 2000-2005*. Clin Infect Dis, 2006. 43(5): p. 585-92.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses

Copyright © 2009 LifeSensors, Inc. All Rights Reserved

- 9) Chiu, S.S., et al., *Human coronavirus NL63 infection and other coronavirus infections in children hospitalized with acute respiratory disease in Hong Kong, China*. Clin Infect Dis, 2005. 40(12): p. 1721-9.
- 10) Bastien, N., et al., *Human coronavirus NL-63 infections in children: a 1-year study*. J Clin Microbiol, 2005. 43(9): p. 4567-73.
- 11) Bastien, N., et al., *Human coronavirus NL63 infection in Canada*. J Infect Dis, 2005. 191(4): p. 503-6.
- 12) Arden, K.E., et al., *New human coronavirus, HCoV-NL63, associated with severe lower respiratory tract disease in Australia*. J Med Virol, 2005. 75(3): p. 455-62.
- 13) Vabret, A., et al., *Human coronavirus NL63, France*. Emerg Infect Dis, 2005. 11(8): p. 1225-9.
- 14) Chen, Z., et al., *Proteolytic processing and deubiquitinating activity of papain-like proteases of human coronavirus NL63*. J Virol, 2007. 81(11): p. 6007-18.
- 15) Nicholson, B., et al., *Characterization of ubiquitin and ubiquitin-like-protein isopeptidase activities*. Protein Sci, 2008.

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