

## Data Sheet

### AC-HLY, His-FLAG-Tags

*B. pertussis*, Recombinant, N-terminal His-FLAG-tag

**Catalog #:** 11005

**Lot #:** 190814-C

**Conc.:** 6.24 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 200 mM Imidazole and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Parkhill, J., *et al.*, *Nature Genetics*, 2003; **35**: 32-40.
2. Althouse, B.M., *et al.*, *BMC Medicine*, 2015 Mar 29; **13**(146).

**Description:** *B. pertussis* Bifunctional Hemolysin/Adenylate Cyclase (AC-HLY), also known as CyaA, Genbank Accession No. NP\_879578, a.a. 2-400 with N-terminal His-FLAG-tag, expressed in an *E. coli* expression system. MW = 46 kDa.

## Quality Assurance

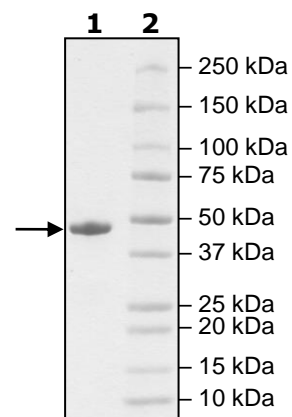
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2  $\mu\text{g}$  AC-HLY

**Lane 2:**  
Protein Marker

**MW:** 46 kDa

**Purity:**  $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)