

## Epitope Tag Small Motif Antibody Sampler Kit

**Catalog# / Size** 699903 / 1 kit

**Regulatory Status** RUO

**Description** The Epitope Tag (Small Motif) Antibody Sampler Kit provides a convenient resource to detect small Epitope Tag fused protein by Western Blot. The DYKDDDDK tag, commonly referred to as Sigma®'s FLAG® Tag, is often used as a protein modification in order to simplify the labeling and detection of proteins. The HA tag (hemagglutinin) is an amino acid sequence derived from the human influenza hemagglutinin surface glycoprotein, corresponding to amino acids 98-106. Anti-His tag antibody was raised against linear 8XHis tag. It reacts with His tag protein, either C, N, or internal His6 and up repeats. The 9E10 monoclonal antibody recognizes human myc and the 10 amino acid epitope tag of human c-myc. The V5 tag is a 14 amino acid peptide derived from a small epitope (Pk) on the P and V proteins of simian virus 5 (SV5), a member of the paramyxovirus family. For additional antigen details, please refer to individual product datasheets.

### Kit Contents

#### Kit Contents

Specificity	Clone	Size	Reactivity	Isotype	MW (kD)
<a href="#">Anti-DYKDDDDK</a>	L5	25 µg	DYKDDDDK tag epitope	Rat IgG2a, κ	N/A
<a href="#">Anti-c-Myc</a>	9E10	25 µg	EQKLISEEDL tag epitope	Mouse IgG1	N/A
<a href="#">Anti-HA.11</a>	16B12	25 µg	YPYDVPDYA tag epitope	Mouse IgG1, κ	N/A
<a href="#">Anti-V5 Tag</a>	7/4	25 µg	GKPIPNPLLGLDST tag epitope	Mouse IgG	N/A
<a href="#">Anti-His Tag</a>	J099B12	25 µg	6x His tag epitope	Mouse IgG1, κ	N/A

\* For detailed information about each specificity, please follow the provided links.

### Product Details

**Formulation** Please refer to individual product datasheets for details.

**Preparation** All antibodies in this kit were purified by affinity chromatography.

**Concentration** All antibodies in the kit are at 0.5mg/ml

**Storage & Handling** Upon receipt, store undiluted at 2-8°C.

**Application** [WB - Quality tested](#)

**Recommended Usage** Each lot of antibodies in this kit is quality control tested by Western Blotting. For Western blotting, the suggested uses of these reagents are as follows:

Anti-DYKDDDDK: 0.1-1 µg/ml  
 Anti-Myc: 0.1-0.5 µg/ml

Anti-HA.11: 0.1-1 µg/ml  
Anti-V5 tag: 0.1-0.5 µg/ml  
Anti-His tag: 0.1-0.2 µg/ml

It is recommended that the reagent be titrated for optimal performance for each application.

## Application References

(PubMed link indicates  
BioLegend citation)

1. Sasaki M, *et al.* 2011. *J. Biol. Chem.* 286:39370. [PubMed](#) (Anti-DYKDDDDK)
2. Sonder SU, *et al.* 2012. *J Immunol.* 188:5906. [PubMed](#) (Anti-DYKDDDDK)
3. Toyo-Oak K, *et al.* 2014. *J Neurosci.* 34:12168. [PubMed](#) (Anti-DYKDDDDK)
4. Das A, *et al.* 2016. *J. Biol. Chem.* 291: 6096 - 6110. [PubMed](#) (Anti-DYKDDDDK)
5. Kondo, S., *et al.* 2011. *J. Virol.* 85:11255. [PubMed](#) (Anti-Myc)
6. Bourgeois-Daigneault MC, *et al.* 2012. *J Immunol.* 188:4959. [PubMed](#) (Anti-Myc)
7. Robu M, *et al.* 2013. *PNAS.* 110:1658. [PubMed](#) (Anti-Myc)
8. Pecot MY, Malhotra V. 2004. *Cell.* 116:99. (Anti-HA.11)
9. Görtz D, *et al.* 2015. *Scientific Reports.* 5: 14685. [PubMed](#) (Anti-HA.11)
10. Thoms M, *et al.* 2016. *Nucleic Acids Res.* 44: 926 - 939. [PubMed](#) (Anti-HA.11)
11. Burritt JB, *et al.* 1995. *J. Biol. Chem.* 270:16974. (Anti-His tag)
12. Taylor RM, *et al.* 2007. *Methods Mol. Biol.* 412:429. (Anti-His tag)
13. Keshwani M, *et al.* 2015. *Biochem. J.* 472: 329 - 338. [PubMed](#) (Anti-His tag)

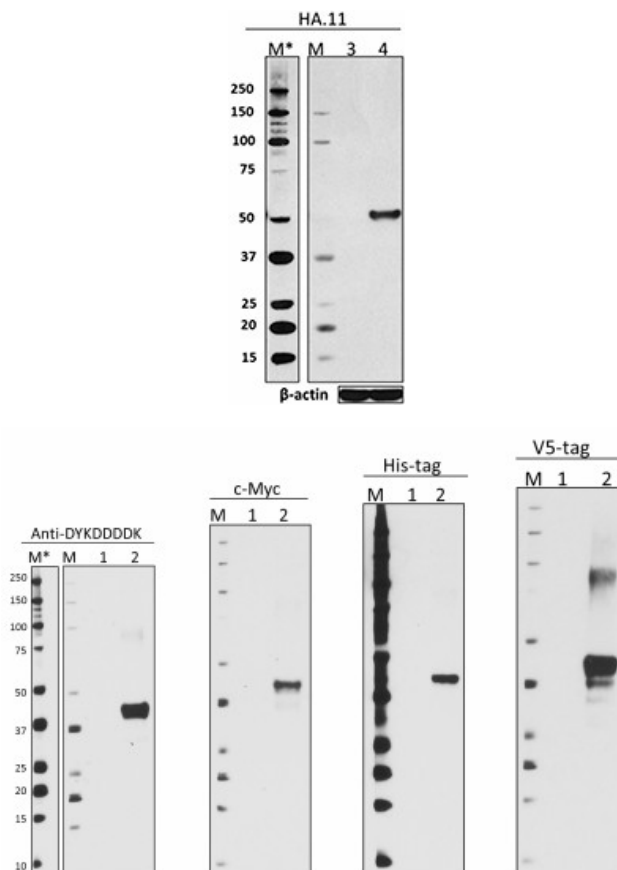
## Antigen Details

**Biology Area** Cell Biology  
**Gene ID** [4609](#)

## Related Protocols

[Western Blotting Protocol](#)

## Product Data



15µg each of total protein from CHO cells (lane 3), and CHO transfected with HA tagged protein (lane 4) were resolved by 4-12% Bis-tris gel electrophoresis, transferred to nitrocellulose, and probed with 1 µg/mL Purified anti-HA.11 Epitope Tag (clone 16B12) Antibody. Proteins were visualized using an HRP goat anti-mouse IgG (clone Poly4053) secondary antibody and chemiluminescence detection. Direct-Blot™ HRP anti-β-actin (clone 2F1-1) Antibody was used as a loading control. Lane M: MW ladder. Lane M\*: long exposure MW ladder.

50ng recombinant human IFN-γ protein (lane 1, negative control) and Posi-Tag Epitope Tag Protein (lane 2) were resolved by 4-12% Bis-tris gel electrophoresis, transferred to nitrocellulose, and probed with either 1 µg/mL of Purified anti-DYKDDDDK Tag (clone L5) Antibody, 0.5 µg/ml of anti-C-Myc (clone 9E10) antibody, 0.2 µg/mL of Purified anti-His Tag (clone J099B12) or anti-V4-tag (clone 7/4) antibody. Proteins were visualized using an HRP goat anti-mouse IgG (clone Poly4053) secondary antibody and chemiluminescence detection. Lane M: MW ladder. Lane M\*: long exposure MW ladder.

For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

\*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, [www.biolegend.com/ordering#license](http://www.biolegend.com/ordering#license)). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

8999 BioLegend Way, San Diego, CA 92121 [www.biolegend.com](http://www.biolegend.com)  
Toll-Free Phone: 1-877-Bio-Legend (246-5343) Phone: (858) 768-5800 Fax: (877) 455-9587