

---

## FLAG Tag Antibody

(Catalog No. A26003)

---

### Description

FLAG Tag Antibody is specific for the hydrophilic epitope tagging octapeptide sequence DYKDDDDK (Asp-Tyr-Lys-Asp-Asp-Asp-Asp-Lys), when fused to either the amino- or carboxyl- terminus or at an internal site on target proteins. Clone 6F7 works at neutral pH and has very low background staining in immunocytochemistry.

### Antibody Type

Monoclonal Antibody

### Purification

Purified IgG - liquid

### Immunogen

### Isotype

IgG1

### Size

100 µg

### Formulation

Phosphate buffered saline. 0.09% Sodium Azide (NaN<sub>3</sub>)

### Specificity

Synthetic Peptide

### Storage

Store at +4°C or at -20°C if preferred. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody.

### Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

### Applications & Suggested Dilutions

ELISA; Flow Cytometry; Immunofluorescence; Immunohistology - Paraffin; Immunoprecipitation; Western Blotting

### Ordering Information:

Products	Size	Cat. No.
FLAG Tag Antibody	100 µg	A26003-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*