

**FITC Mouse Anti-human/mouse VCP
Antibody *Hs-14, monoclonal***Catalog number: V1032425
Unit size: 50 tests**Product Details**

| | |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | Lot specific (please consult certificate of analysis for given lot) |
| Formulation | Stabilizing Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA |

Antibody Properties

| | |
|--------------------|--------------|
| Species Reactivity | Human, mouse |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Mouse |
| Immunogen | VCP |
| Clone | Hs-14 |
| Conjugate | FITC |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions |
| Application | FC (QC TESTED) |

Applications

This is an antibody that targets VCP.