

**Purified Rabbit Anti-species independent  
Dendra2 Antibody \*PAb (836), monoclonal\***Catalog number: V103575  
Unit size: 0.1 mg**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        | Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

|                    |                     |
|--------------------|---------------------|
| Species Reactivity | Species independent |
| Class              | Primary             |
| Clonality          | Monoclonal          |
| Host               | Rabbit              |
| Immunogen          | Dendra2             |
| Clone              | PAb (836)           |

**Biological Properties**

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

**Applications**

Identifies the photoswitchable Dendra2 tag, a protein derived from Dendronephthya.