

**Purified Mouse Anti-human AGPS Antibody**  
**\*AGPS-03, monoclonal\***Catalog number: V103040  
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	AGPS
Clone	AGPS-03

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC, WB

**Applications**

---

Alkyldihydroxyacetonephosphate synthase, peroxisomal (sometimes called aging-associated gene 5 protein) is a protein with a molecular weight of 85 kDa, located in the peroxisomal membrane, peroxisome and cytosol of cells. Sequencing of alkyldihydroxyacetonephosphate synthase, peroxisomal has shown it contains a primary structural unit, the FAD-binding PCMH-type domain. Alkyldihydroxyacetonephosphate synthase, peroxisomal has been thought to be involved with essential functions like alkyglycerone-phosphate synthase and oxidoreductase activity, and it recognizes FAD. Alkyldihydroxyacetonephosphate synthase, peroxisomal aids in organismal processes, for instance, ether lipid biosynthetic process. Alkyldihydroxyacetonephosphate synthase, peroxisomal is clinically significant because abnormalities in its function have been closely linked to diseases like rhizomelic chondrodysplasia punctata 3 (RCDP3). Rhizomelic chondrodysplasia punctata 3, an autosomal recessive inheritancedisorder characterized by short humerus, epiphyseal stippling and failure to thrive, has in particular been of interest to researchers.