



02-6100 1 mg  
Purified Mouse  
Myeloma IgG<sub>1</sub>  
Lot No. See product label

### MOUSE MYELOMA IgG<sub>1</sub>

**FORM:**

Liquid: 1 ml.

**COMPOSITION:**

This purified IgG<sub>1</sub> protein is in 10 mM phosphate-buffered saline, pH 7.4, containing 0.1% sodium azide as a preservative. The protein concentration is 1 mg/ml based on E<sub>(1%)</sub> =14 at 280 nm.

**SOURCE:**

This protein was purified from mouse myeloma IgG<sub>1</sub> ascites induced by myeloma cells.

**PURITY:**

This mouse IgG<sub>1</sub> was characterized by ELISA.

**STORAGE:**

Store at 2-8°C for up to one month. Store at -20°C for long term storage. Avoid repeated freezing and thawing.

**For Research Use Only**

SS001117

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288  
E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI026100

(Rev 12/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). 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.



02-6100 1 mg  
Purified Mouse  
Myeloma IgG<sub>1</sub>  
Lot No. See product label

### MOUSE MYELOMA IgG<sub>1</sub>

**FORM:**

Liquid: 1 ml.

**COMPOSITION:**

This purified IgG<sub>1</sub> protein is in 10 mM phosphate-buffered saline, pH 7.4, containing 0.1% sodium azide as a preservative. The protein concentration is 1 mg/ml based on E<sub>(1%)</sub> =14 at 280 nm.

**SOURCE:**

This protein was purified from mouse myeloma IgG<sub>1</sub> ascites induced by myeloma cells.

**PURITY:**

This mouse IgG<sub>1</sub> was characterized by ELISA.

**STORAGE:**

Store at 2-8°C for up to one month. Store at -20°C for long term storage. Avoid repeated freezing and thawing.

**For Research Use Only**

SS001117

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288  
E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI026100

(Rev 12/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). 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.