



33-5700 100 µg
 Mouse Anti-Osteocalcin
 Lot No.

MOUSE MONOCLONAL ANTI-OSTEOCALCIN

FORM

Lyophilized. This antibody was purified by ammonium sulfate fractionation and ion-exchange chromatography from mouse ascites fluid, dissolved in 0.15 M PBS (pH 7.4) containing 1.0% bovine serum albumin and 0.1% sodium azide. Reconstitute with 50 µl distilled water to give a 2 mg/ml stock solution.

CLONE: OCG-3

ISOTYPE: IgG₃

IMMUNOGEN

Bovine Osteocalcin⁽¹⁾. Monoclonal antibody was obtained by fusing the mouse myeloma cell-line P3U1 with spleen cells of BALB/c mouse after immunization with bovine Osteocalcin⁽¹⁾.

SPECIFICITY

OCG-3 recognizes γ-carboxylated residues 21-31 (does not react with heat decarboxylated Osteocalcin).

REACTIVITY

This product reacts with both bovine and human Osteocalcin.

SAMPLE	ELISA	WB
Bovine	+	+
Human	+	+

BACKGROUND

This antibody can be used effectively for research of metabolic bone disease and ectopic calcification.

USAGE

The dilutions listed below are good starting points; however, optimal dilution of the antibody should be determined by the investigator for each application.

ELISA: 1:1000 (2 µg/ml)

Western blotting: 2-10 µg/ml

(cont'd)

www.invitrogen.com

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

PI335700

(Rev 10/08) 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). 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.



33-5700 100 µg
 Mouse Anti-Osteocalcin
 Lot No.

MOUSE MONOCLONAL ANTI-OSTEOCALCIN

FORM

Lyophilized. This antibody was purified by ammonium sulfate fractionation and ion-exchange chromatography from mouse ascites fluid, dissolved in 0.15 M PBS (pH 7.4) containing 1.0% bovine serum albumin and 0.1% sodium azide. Reconstitute with 50 µl distilled water to give a 2 mg/ml stock solution.

CLONE: OCG-3

ISOTYPE: IgG₃

IMMUNOGEN

Bovine Osteocalcin⁽¹⁾. Monoclonal antibody was obtained by fusing the mouse myeloma cell-line P3U1 with spleen cells of BALB/c mouse after immunization with bovine Osteocalcin⁽¹⁾.

SPECIFICITY

OCG-3 recognizes γ-carboxylated residues 21-31 (does not react with heat decarboxylated Osteocalcin).

REACTIVITY

This product reacts with both bovine and human Osteocalcin.

SAMPLE	ELISA	WB
Bovine	+	+
Human	+	+

BACKGROUND

This antibody can be used effectively for research of metabolic bone disease and ectopic calcification.

USAGE

The dilutions listed below are good starting points; however, optimal dilution of the antibody should be determined by the investigator for each application.

ELISA: 1:1000 (2 µg/ml)

Western blotting: 2-10 µg/ml

(cont'd)

www.invitrogen.com

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

PI335700

(Rev 10/08) 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). 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.

(33-5700 cont'd)

RELATED PRODUCTS:

Goat x Mouse IgG (H+L)

Conjugate	Cat. No.
Purified	62-6500
FITC	62-6511
TRITC	62-6514
Cy TM 3	62-6515
Cy TM 5	62-6516
HRP	62-6520
Alk. Phos.	62-6522
Biotin	62-6540
Sepharose [®] -4B	62-6541
Goat x Mouse IgGAM (H+L) (DS grade)	Purified 65-6400
	FITC 65-6411
	HRP 65-6420
	Alk. Phos. 65-6422
	Biotin 65-6440

STORAGE

Store at -20°C for long term storage. Store in aliquots, avoid repeated freezing and thawing.

REFERENCES

1. Hauschka, P.U., Lian, J.B., Gallop, P.M., (1975) *Proc. Natl. Acad. Sci. U.S.A.*,72:3925.

NOT FOR DIAGNOSTIC USE

(33-5700 cont'd)

RELATED PRODUCTS:

Goat x Mouse IgG (H+L)

Conjugate	Cat. No.
Purified	62-6500
FITC	62-6511
TRITC	62-6514
Cy TM 3	62-6515
Cy TM 5	62-6516
HRP	62-6520
Alk. Phos.	62-6522
Biotin	62-6540
Sepharose [®] -4B	62-6541
Goat x Mouse IgGAM (H+L) (DS grade)	Purified 65-6400
	FITC 65-6411
	HRP 65-6420
	Alk. Phos. 65-6422
	Biotin 65-6440

STORAGE

Store at -20°C for long term storage. Store in aliquots, avoid repeated freezing and thawing.

REFERENCES

2. Hauschka, P.U., Lian, J.B., Gallop, P.M., (1975) *Proc. Natl. Acad. Sci. U.S.A.*,72:3925.

NOT FOR DIAGNOSTIC USE

www.invitrogen.com

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

PI335700

(Rev 10/08) 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). 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.

www.invitrogen.com

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

PI335700

(Rev 10/08) 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). 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.