

Collagen I Matrix Solution (rat)

Collagen I Matrix for 3D cell culture

Instruction Manual

Catalog Number	PK-CA577-K916-5
Description	<p>Collagen is the most widely used extracellular matrix (ECM) protein for cell culture. It is involved in cell attachment, growth/proliferation, differentiation, migration, and tissue morphogenesis during development.</p> <p>Collagen I is the most common fibrillar collagen and is the major structural component of extracellular matrices found in connective tissue and internal organs.</p> <p>PromoCell's high-quality Collagen I is purified from rat tail tendons. It is capable to polymerize, forming a hydrogel at 37°C and directing cells to grow in a 3D fashion and assemble into organotypic structures.</p> <p>PromoCell's Collagen I provides in vivo-like extracellular matrix (ECM) structures. It is usable for the preparation of a variety of gel concentrations and shows a fast polymerization that facilitates optimal cell distribution in 3D gels.</p> <p>It is a component of our Cell Invasion Kits (Cat. No. PK-CA577-K916 & PK-CA577-K917).</p>
Quantity	5 ml
Applications	<ul style="list-style-type: none">• 3D cell culture• Cell invasion and angiogenesis assays• Cell adhesion studies
Applications / Assay Protocol	For Cell Invasion Assay please see protocol of kits PK-CA577-K916 & PK-CA577-K917.
Intended Use	For in vitro research use only. Not for diagnostic or therapeutic procedures.
Storage & Stability	Store at -20°C, protected from light. Briefly centrifuge vial prior to opening. Aliquot under hood and store at 4°C, if needed.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

USA/Canada

Phone: 1 - 866 - 251 - 2860 (toll free)
Fax: 1 - 866 - 827 - 9219 (toll free)

Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)
Fax: 0800 - 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 - 96 03 33 (toll free)
Fax: 0800 - 169 85 54 (toll free)

Other Countries

Phone: +49 6221 - 649 34 0
Fax: +49 6221 - 649 34 40