

TECHNICAL MANUAL

Unpacking and Handling

Catalogue number: MAB-802-VIAL/ 1 vial of ES-Collect™
Y20011

1. Unpack ES-Collect™ as soon as possible after arrival.
2. Check all materials for leakage or breakage.
3. ES-Collect™ should be stored at +2-8°C upon arrival.

For technical support email: tech-cellartis@takara-clontech.eu

Authorised uses

Except as otherwise agreed in writing, the purchase of goods only conveys to you the non-transferable right for only you to use the quantity of goods and components of goods purchased in compliance with the applicable intended use statement. Unless otherwise authorized, no right to resell the goods, or any portion of them, is conveyed hereunder.

The goods are intended for research use only and are not to be used for any other purposes including, but not limited to: unauthorized commercial purposes, *in vitro* diagnostic purposes, *ex vivo* or *in vivo* therapeutic purposes, investigational use, in foods, drugs, devices or cosmetics of any kind, or for consumption by or use in connection with or administration or application to humans or animals.