

ab286892 – Lymphoprep

For the isolation of human mononuclear cells.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286892>

Storage and Stability

Store at +4°C to +30°C. Protect from air and light. Prolonged exposure to direct sunlight leads to release of iodine from the sodium diatrizoate molecule.

Protocol

1. Collect blood into a tube containing anticoagulant (EDTA, heparin, ACD) or use defibrinated blood.
2. Dilute the blood by adding an equal volume of 0.9% NaCl.
3. Carefully layer 6 ml of the diluted blood over 3 ml Lymphoprep in a 12-15 mm centrifuge tube. Alternatively, Lymphoprep can be underlayered. Avoid mixing of blood and separation fluid. Cap the tube to prevent the formation of aerosols.
4. Centrifuge at 800 x g for 20 min at RT (approximately 20°C) in a swing-out rotor. If the blood is stored for more than 2 hr increase the centrifugation time to 30 min.
5. After centrifugation the mononuclear cells form a distinct band at the sample/medium interface, as shown in the figure. The cells are best removed from the interface using a Pasteur pipette without removing the upper layer.
6. Dilute the harvested fraction with 0.9% NaCl or medium to reduce the density of the solution and pellet the cells by centrifugation for 10 min at 250 x g.

Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)

Version 1, Last updated Friday, October 15, 2021