

Version 1: Last updated 19 February 2018

ab105229

Discovery XT[®] Staining Protocol for J-chain on Human Tonsil

Clone ID: SP105 Host Species: Rabbit Monoclonal

This product is for research use only and is not intended for diagnostic use.

Procedure: Res IHC Omni-UltraMap HRP XT

1. Select Tissue Sample
2. Select Paraffin
3. Select Deparaffinization
4. Select Cell Conditioning
5. Select Conditioner #1
6. Select Mild CC1
7. Select Antibody
8. Select Antibody Titration
9. Select Standard Titration Ab Inc
 - Instrument will hold before Antibody Titration, press the button on the instrument when prompted on screen to continue to Antibody Titration step
 - Open slide drawer and manually apply 100ul of antibody diluted 1:50 in Antibody Diluent (ab227819)
 - Close slide drawer and press the button on the instrument to continue run
10. Incubate Primary Antibody for [32 Minutes] [with Heat: default setting]
11. Select [OmniMap anti-Rb HRP] (Multimer HRP) and Incubate for [8 Minutes]
12. Select Counterstain
13. Select Standard
 - Instrument will hold before Counterstain, press the button on the instrument when prompted on screen to continue to Counterstain step

14. Select [HEMATOXYLIN II] (Counterstain) and Incubate for [4 Minutes]

15. Select Post Counterstain

16. Select [BLUING REAGENT] (Post Counterstain), Incubate for [4 Minutes]

Note: This is a suggested starting point for use of this antibody on DXT staining platform, further optimization may be required.

Technical Support

All information / detail is correct at time of going to print. Technical support questions, please contact Abcam Representatives.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | +86-21-5110-5938 | 400-628-6880

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abcam.com | +64-(0)9-909-7829