

Version 1: Last updated 19 February 2018

ab228462

Discovery XT[®] Staining Protocol for
PD-L1/CD274 on Human Non-Small
Cell Lung Cancer

Clone ID: SP142 Host Species: Rabbit Monoclonal

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

Procedure: Res Hapten IHC XT v1

1. Select Tissue Sample
2. Select Paraffin
3. Select Deparaffinization
4. Select Cell Conditioning (CC1 56 minutes)
5. Select Inhibitor
6. Select Antibody
7. Select antibody titration
8. 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 1:25 in Antibody Diluent (ab227819)
 - Close slide drawer and press the button on the instrument to continue run
9. Select Disable heat for 1st Antibody
10. Select [PREP KIT 1](Antibody) and incubate for [32 mins] (1:25 dilution)
11. Select Linking Antibody
12. Select 2nd Antibody
13. Warmup Slide to [36 Deg C] from Very Low Temperatures (2nd Antibody)

14. Select [Anti-Rabbit HQ] (Detection #1) and incubate for [12Min]
15. Select Enzyme conjugate
16. Select [Anti-HQ HRP] (Conjugate #1), and Incubate for [12 Minutes]
17. Select DAB
18. Select Counterstain
19. Select Use RB for Counterstain
20. Select [HEMATOXYLIN II] (Counterstain) and Incubate for [4 Mins]
21. Select Post Counterstain
22. Select Use RB for Post Counterstain
23. 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