

Anti-Drug Antibody (ADA) Bridging ELISA

ADA - Panitumumab

For use with anti-panitumumab monoclonal antibody products HCA265 or HCA266

This method provides a procedure for generating an ADA ELISA standard curve with anti-panitumumab antibodies, product codes HCA265 or HCA266. The method should always be used in conjunction with the product and batch specific information provided with each vial (see product datasheets). This protocol will need to be adjusted for use with different detection methods and immunoassay technology platforms.

Reagents

- BSA
- HISPEC immunoassay diluent (BUF049)
- Human Serum (Sigma-Aldrich, H4522)
- Lynx Rapid HRP Antibody Conjugation Kit® (LNK001P-LNK006P)
- PBS
 - 136 mM NaCl
 - 2.68 mM KCl
 - 8.1 mM Na₂HPO₄
 - 1.46 mM KH₂PO₄
- PBST
 - PBS with 0.05% Tween®-20
- QuantaBlu™ fluorogenic peroxidase substrate (Thermo Fisher Scientific, 15169)

Materials

- 384-well microtiter plate, black, square flat-bottom wells, MaxiSorp™ PS (Thermo Fisher Scientific, 460518)
- Fluorescence plate reader

96-well plates can be used instead of 384-well plates, e.g. black, flat-bottom MaxiSorp PS (Thermo Fisher Scientific, 437111). For the 96-well format, use 100 µl (instead of 20 µl) of antigen, antibodies or substrate, and 300 µl for the blocking step.

Method

1. Prepare detection antibody: conjugate panitumumab antibody using a Lynx Rapid HRP Antibody Conjugation Kit.

2. Prepare the panitumumab at 1 µg/ml in PBS. Coat the required number of wells of a 384-well microtiter plate with 20 µl per well of the prepared panitumumab, and incubate overnight at 4°C.
3. Wash the microtiter plate five times with PBST.
4. Block the microtiter plate by adding 100 µl 5% BSA in PBST to each well, and then incubate for 1 hour at room temperature (RT).
5. Wash the microtiter plate five times with PBST.
6. For the standard curve, prepare a dilution series of the anti-panitumumab antibody HCA265 (clone AbD23897_hlgG1) or HCA266 (clone AbD23895_hlgG1) in 10% human serum in PBST in triplicate. Final concentrations of anti-panitumumab antibody should cover the range from 0.1 ng/ml to 10,000 ng/ml. Include a zero anti-panitumumab concentration as the background value.
7. Add 20 µl of anti-panitumumab antibody dilution per well (in triplicate for each standard recommended) and incubate for 1 hour at RT.
8. Wash the microtiter plate five times with PBST.
9. To each well add 20 µl HRP conjugated panitumumab diluted to 2 µg/ml in HISPEC buffer and incubate for 1 hour at RT.
10. Wash the microtiter plate ten times with PBST.
11. Add 20 µl QuantaBlu to each well and measure the fluorescence after 30 minutes.

MaxiSorp™ and QuantaBlu™ are trademarks of Thermo Fisher Scientific
Tween® is a registered trademark of ICI Americas Inc.

V1.11.2015