



Staining solution for specific one-step detection of His-tag fusion proteins!

The His-Tag Detect solution is a ready-to-use reagent for high speed detection of his-tagged proteins.

Application:

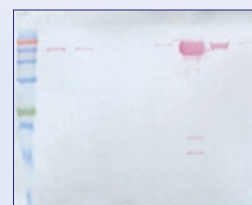
- Detection of His-tagged proteins from samples in SDS-gel:



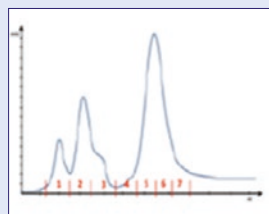
Gel transfer on membrane



+ His-Detect



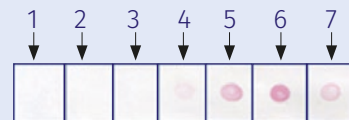
- Detection of fractions containing His-tagged proteins:



Spot fractions on membrane



+ His-Detect



- Expression level analysis of His-tagged proteins at the colony level:



Colony filtration blot1



+ His-Detect



Product

His Detect Solution

Cat.-No.

PS-110

Amount

up to 3 x 3 reactions (30 ml)

