

## Dako Platform Protocol Summary

### IHC Reagents (Dako Platform):

- Envision FLEX Target Retrieval Solution, High pH (Dako, Cat# DM828)
- EnVision FLEX Wash Buffer (Dako, Cat# K8007)
- Envision FLEX Peroxidase-Blocking Reagent (Dako, Cat# SM801)
- Envision FLEX Antibody Diluent (Dako, Cat# S0809)
- Envision FLEX HRP (Dako, Cat# SM802)
- Envision FLEX+ Mouse (Linker) (Dako, Cat# SM804)
- Envision Flex Substrate Buffer (Dako, Cat# SM803)
- Envision Flex DAB+ Chromogen (Dako, Cat# DM827)
- Envision FLEX Hematoxylin (Dako, Cat# SM806)

### Protocol Summary:

#### FLEX+ Mouse

	Category	Code	Reagent Name	Incubation
1	Rinse		Buffer	
2	Enzyme Pretreatment			
3	Rinse		Buffer	
4	Endogenous Enzyme Block	SM801	FLEX Peroxidase Block	5
5	Rinse		Buffer	
6	Primary Antibody			
7	Rinse		Buffer	
8	Secondary Reagent	SM804	FLEX+ Mouse (LINKER)	15
9	Rinse		Buffer	
10	Labelled Polymer	SM802	FLEX /HRP	20
11	Rinse		Buffer	
12	Rinse		Buffer	5
13	Substrate-Chromogen	SM803	FLEX DAB+ Sub-Chromo	10
14	Rinse		Buffer	
15	Counterstain	SM806	FLEX Hematoxylin	5
16	Rinse		DI Water	
17	Rinse		Buffer	5
18	Rinse		DI Water	