



EnVision™ FLEX (Link)
Optional Reagents for Autostainer Link Instruments

Code K8004, K8005, K8006, K8007, K8008, K8009, K8021

These reagents are described in the package inserts for:

Code K8000: EnVision™ FLEX, High pH ([Link](#))

Code K8002 : EnVision™ FLEX+, High pH ([Link](#))

Code K8023: EnVision™ FLEX Mini kit, High pH ([Link](#))

EnVision™ FLEX (Dako Autostainer/Autostainer Plus)
Optional Reagents for Dako Autostainer Instruments

Code K8004, K8005, K8006, K8007, K8018, K8019, K8022

These reagents are described in the package inserts for:

Code K8010: EnVision™ FLEX, High pH (Dako Autostainer/Autostainer Plus)

Code K8012: EnVision™ FLEX+, High pH (Dako Autostainer/Autostainer Plus)

Code K8024: EnVision™ FLEX Mini kit, High pH (Dako Autostainer/Autostainer Plus)