

AbFlex Antibodies Application Data

The table below references applications for which individual AbFlex recombinant antibodies have been validated. To learn more about AbFlex antibodies and our Sortag labeling kits, visit us online at www.activemotif.com/abflex

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	61479	91187	5mC
MeDIP	+	NT	
DB	+	+	
	39937	91217	Ago 1/2/3
ChIP	+	NT	
WB	+	NT	
IP	+	+	
IF	+	NT	
IHC	+	NT	
ICC	+	NT	
	61751	91111	AM Tag
WB	+	+	
ChIP	+	+	
ELISA	NT	+	
	61751	91109	AM Tag (biotin)
ChIP	+	+	
WB	+	+	
ELISA	NT	+	
	39529	91207	ATM phospho Ser1981
WB	+	+	
DB	NT	+	
IF	+	NT	
ICC	+	NT	
	39993	91195	Bmi-1
ChIP	+	NT	
WB	+	+	
IF	+	NT	
IHC	+	NT	
	61577	91123	Cas9
WB	+	+	
IF	+	NT	

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	61577	91123	Cas9
ICC	+	NT	
IP	+	NT	
ELISA	NT	+	
	61203	91135	EED
ChIP	+	+	
ChIP-Seq	+	+	
IHC	+	NT	
Bead ELISA	NT	+	
	39531	91215	HDAC1
ChIP	+	NT	
WB	+	+	
IP	+	NT	
IF	+	NT	
IHC	+	NT	
ICC	+	NT	
	39533	91197	HDAC2
ChIP	+	+	
ChIP-Seq	NT	+	
WB	+	+	
IF	+	NT	
ICC	+	NT	
	61275	91139	Hif-1 alpha
WB	+	NT	
IHC	+	NT	
ELISA	NT	+	
	61251	91103	H3K9ac
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IF	+	+	
DB	+	NT	
ICC	+	+	
ELISA	NT	+	

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	61251	91101	H3K9ac (biotin)
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IF	+	+	
DB	+	NT	
ICC	+	+	
ELISA	NT	+	
	61061	91127	H3K56ac
ChIP	+	+	
ChIP-Seq	NT	+	
DB	+	NT	
ELISA	NT	+	
	39671	91107	H4K20me3
ChIP	+	+	
ChIP-Seq	NT	+	
WB	+	+	
IF	NT	+	
ICC	NT	+	
DB	+	NT	
ELISA	NT	+	
	39671	91105	H4K20me3 (biotin)
ChIP	+	+	
ChIP-Seq	NT	+	
WB	+	+	
IF	NT	+	
ICC	NT	+	
DB	+	NT	
ELISA	NT	+	
	39815	91143	JMJD2A
WB	+	+	
ELISA	NT	+	
	39287	91189	Lamin A/C
WB	+	+	
IF	+	+	
ICC	+	NT	

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	39935	91205	MAZ
ChIP	+	NT	
WB	+	+	
EMSA	+	NT	
	39789	91201	MITF
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	NT	
IP	+	NT	
IF	+	NT	
EMSA	+	NT	
ICA	+	NT	
	39279	91203	Myc-Tag
WB	+	+	
	61075	91213	c-Myc
WB	+	+	
IP	+	NT	
EMSA	+	NT	
	61147	91243	Notch1
WB	+	+	
	39541	91213	Nucleolin
WB	+	+	
IF	+	NT	
ICC	+	NT	
	39553	91247	p53
ChIP	+	NT	
WB	+	+	
IP	+	NT	
	61155	91227	PBX2
WB	+	+	
IHC	+	NT	
	39383	91245	Rad21
WB	+	+	
IF	+	+	
ICC	+	NT	

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	39097	91151	RNA Pol II
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IF	+	+	
ICC	+	NT	
	61383	91219	RNA Pol II CTD phospho Tyr1
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IP	+	NT	
IF	+	+	
ELISA	NT	+	
ICC	+	+	
	61083	91115	RNA Pol II CTD phospho Ser2
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IP	+	NT	
IF	+	+	
ELISA	NT	+	
ICC	+	+	
	61085	91119	RNA Pol II CTD phospho Ser5
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IP	+	NT	
IF	+	+	
ELISA	NT	+	
ICC	+	+	
	61085	91117	RNA Pol II CTD phospho Ser5 (biotin)
ChIP	+	+	
ChIP-Seq	+	+	
WB	+	+	
IP	+	NT	
IF	+	+	

Application	Parent Antibody	AbFlex Antibody	Antibody Target
	61085	91117	RNA Pol II CTD phospho Ser5 (biotin)
ELISA	NT	+	
ICC	+	+	
	61465	91225	SMARCA1 / SNF2L1
WB	+	+	
IHC	+	NT	
	61741	91171	TET1
ChIP	+	NT	
ChIP-Seq	+	NT	
WB	+	+	
IP	+	NT	
IF	+	NT	
ICC	+	NT	