

アクティブ・モティフ AbFlex®抗体プレゼント！

- 抗体のご購入でサンプルサイズを無料提供します -



キャンペーン期間

～ 2022/6/30

抗体のロット間格差や標識に
悩まされていませんか？

AbFlex®抗体
お試しください！！

AbFlex®抗体の特長

- ・ **ChIP, ChIP-Seq**検証済み抗体多数！
- ・ 様々な標識・精製法に柔軟に対応
Sortase認識モチーフ (LPXTG)・ビオチン化タグ・6xHisタグ
- ・ ロット間で一貫した再現性、信頼性の高い長期的な供給
- ・ 高い特異性と費用対効果
- ・ アニマルフリー生産



AbFlex®抗体
詳細はコチラ

モノクローナル抗体、ポリクローナル抗体 (100 µg または 100 µlサイズ) ご注文の際に、ご希望の AbFlex®リコンビナント抗体 (サンプルサイズ/10 µg) のカタログ番号をお知らせいただきますと、無料でご提供します。

その他の割引との併用はできません。また、キャンペーン開始前のご購入には適用されません。抗体1個のご注文につきサンプル1個とさせていただきます。

mAb, pAb (100µg scale) のご注文
ご希望のAbFlex®抗体 (Cat#) を併記

AbFlex®抗体 プレゼント！
(10µg)

プレゼント対象 AbFlex® 抗体は裏面へ

プレゼント対象 AbFlex®リコンビナント抗体 (サンプルサイズ 10 µg)

製品名	Cat#	製品名	Cat#
AbFlex 5-methylcytosine antibody (rAb)	91188	AbFlex HSP72 antibody (rAb)	91304
AbFlex Ago 1/2/3 antibody (rAb)	91218	AbFlex ING5 antibody (rAb)	91330
AbFlex AM-Tag antibody (rAb) *	91112	AbFlex JMJD2A antibody (rAb)	91144
AbFlex ATM phospho Ser1981 antibody (rAb)	91208	AbFlex JMJD2B / KDM4B antibody (rAb)	91306
AbFlex Bcl6 antibody (rAb)	91288	AbFlex JMJD2C / KDM4C antibody (rAb)	91308
AbFlex BMI-1 antibody (rAb) *	91196	AbFlex JMJD2D / KDM4D antibody (rAb)	91370
AbFlex BRD4 antibody (rAb) *	91302	AbFlex KDM1B / LSD2 antibody (rAb)	91328
AbFlex BrdU antibody (rAb)	91270	AbFlex KDM5A antibody (rAb) *	91212
AbFlex c-Myc antibody (rAb)	91214	AbFlex Lamin A/C antibody (rAb)	91190
AbFlex Cas9 antibody (rAb) *	91124	AbFlex MAZ antibody (rAb)	91206
AbFlex CHD3 antibody (rAb)	91274	AbFlex MEIS 1/2/3 antibody (rAb)	91260
AbFlex CTCF antibody (rAb) *	91286	AbFlex METTL3 antibody (rAb)	91334
AbFlex Dicer antibody (rAb)	91242	AbFlex Mi-2 beta (CHD4) antibody (rAb)	91276
AbFlex DNMT1 antibody (rAb)	91176	AbFlex Mili / PiwiL2 antibody (rAb)	91254
AbFlex EED antibody (rAb) *	91136	AbFlex MITF antibody (rAb) *	91202
AbFlex EHMT2 (G9a) antibody (rAb)	91230	AbFlex MPP8 antibody (rAb)	91278
AbFlex EZH2 antibody (rAb)	91160	AbFlex Myc-Tag antibody (rAb)	91204
AbFlex HA-Tag antibody (rAb)	91272	AbFlex N6-Methyladenosine (m6A) antibody (rAb)	91262
AbFlex HDAC1 antibody (rAb) *	91216	AbFlex Nanog antibody (rAb)	91252
AbFlex HDAC2 antibody (rAb) *	91198	AbFlex Notch1 antibody (rAb)	91244
AbFlex HDAC3 antibody (rAb)	91200	AbFlex Nucleolin antibody (rAb)	91232
AbFlex HIF-1 alpha antibody (rAb)	91140	AbFlex p53 antibody (rAb)	91248
AbFlex His-Tag antibody (rAb)	91184	AbFlex PBX2 antibody (rAb)	91228
AbFlex Histone H2A antibody (rAb)	91326	AbFlex Pim2 antibody (rAb)	91240
AbFlex Histone H2A.XS139ph antibody (rAb)	91396	AbFlex Rad21 antibody (rAb)	91246
AbFlex Histone H3, C-terminal antibody (rAb, Mouse)	91298	AbFlex RNA Pol II antibody (rAb) *	91152
AbFlex Histone H3, C-terminal antibody (rAb, Rabbit) *	91300	AbFlex RNA pol II CTD phospho Ser2 antibody (rAb) *	91116
AbFlex Histone H3.3 antibody (rAb) *	91192	AbFlex RNA pol II CTD phospho Ser5 antibody (rAb) *	91120
AbFlex Histone H3K14ac antibody (rAb)	91398	AbFlex RNA Pol II CTD phospho Tyr1 antibody (rAb) *	91220
AbFlex Histone H3K27ac antibody (rAb) *	91194	AbFlex SaCas9 antibody (rAb)	91280
AbFlex Histone H3K27me3 antibody (rAb)	91168	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM001414)	91376
AbFlex Histone H3K27me3 antibody (rAb) *	91404	(Clone AM001414 (414-1))	
AbFlex Histone H3K36me3 antibody (rAb) *	91266	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM002414)	91364
AbFlex Histone H3K4me0 antibody (rAb, rabbit)	91318	(Clone AM002414 (414-2))	
AbFlex Histone H3K4me0 antibody (rAb)	91268	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM004414)	91374
AbFlex Histone H3K4me1 antibody (rAb)	91290	(Clone AM004414 (414-4))	
AbFlex Histone H3K4me2 antibody (rAb) *	91322	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM009105)	91366
AbFlex Histone H3K4me3 antibody (rAb) *	91264	(Clone AM009105 (105-9))	
AbFlex Histone H3K56ac antibody (rAb) *	91128	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM038105)	91372
AbFlex Histone H3K79me2 antibody (rAb) *	91400	(Clone AM038105 (105-38))	
AbFlex Histone H3K9ac antibody (rAb) *	91104	AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM043105)	91368
AbFlex Histone H3K9me0 antibody (rAb) *	91156	(Clone AM043105 (105-43))	
AbFlex Histone H3K9me2 antibody (rAb)	91336	AbFlex SMARCA1 / SNF2L1 antibody (rAb)	91226
AbFlex Histone H3S10ph antibody (rAb)	91132	AbFlex Snail antibody (rAb)	91256
AbFlex Histone H4 antibody (rAb)	91296	AbFlex Supt5H/SPT5 antibody (rAb)	91332
AbFlex Histone H4K20me1 antibody (rAb)	91292	AbFlex Suz12 antibody (rAb)	91164
AbFlex Histone H4K20me3 antibody (rAb) *	91108	AbFlex TET1 antibody (rAb) *	91172
AbFlex Histone H4K5ac/K8ac antibody (rAb) *	91320	AbFlex TWIST antibody (rAb) *	91324
		AbFlex Ubiquitin antibody (rAb)	91258

* ChIP or ChIP-Seq 検証済み



アクティブ・モティフ株式会社

〒162-0824 東京都新宿区揚場町2-21
 Tel: 03-5225-3638 Fax: 03-5261-8733
 E-mail: japantech@activemotif.com
 www.activemotif.com