

一次抗体キャンペーン

2本以下購入 **20% off** 3本以上購入(組み合わせ自由) **25% off**

キャンペーン期間: 2020年1月27日(月)~2020年3月31日(火) 受注分まで

エピジェネティクス研究に重点をおいたアクティブ・モティフ社の一次抗体ラインナップは、ヒストンおよびDNA修飾をはじめ、転写因子、クロマチン修飾因子および幹細胞制御因子まで多岐にわたります。今回のキャンペーンは一次抗体ならば包装や種類(ポリクローナル、モノクローナル、リコンビナント抗体)を問わず、全製品が対象です。

以下に人気の高い製品を掲載します。掲載されていない抗体も多数ございますので、裏面に掲載のURLまたはQRコードをご参照の上、弊社ウェブサイトにてご確認ください。

ヒストン/ヒストン修飾 (200種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
Histone H2A.Z antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	100 µl	39113	¥70,000	→ ¥56,000	¥52,500
Histone H3K27ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, IHC, WB	100 µg	39133	¥70,000	→ ¥56,000	¥52,500
Histone H3K9ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	100 µl	39137	¥70,000	→ ¥56,000	¥52,500
Histone H3ac (pan-acetyl) antibody (pAb)	ChIP, ChIP-Seq, DB, WB	100 µl	39139	¥70,000	→ ¥56,000	¥52,500
Histone H3K4me2 antibody (pAb)	ChIP, ChIP-Seq, DB, WB	100 µl	39141	¥70,000	→ ¥56,000	¥52,500
Histone H3K79me2 antibody (pAb)	ChIP, ChIP-Seq, DB, WB	100 µl	39143	¥70,000	→ ¥56,000	¥52,500
Histone H3K27me3 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, IHC, WB	100 µg	39155	¥70,000	→ ¥56,000	¥52,500
Histone H3K4me3 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, IHC, WB	100 µl	39159	¥70,000	→ ¥56,000	¥52,500
Histone H3K9me3 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, IHC, WB	100 µl	39161	¥70,000	→ ¥56,000	¥52,500
Histone H4K12ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µl	39165	¥70,000	→ ¥56,000	¥52,500
Histone H3K9me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	100 µl	39239	¥70,000	→ ¥56,000	¥52,500
Histone H3K27me2 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	200 µl	39245	¥70,000	→ ¥56,000	¥52,500
Histone H3K36me2 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µl	39255	¥70,000	→ ¥56,000	¥52,500
Histone H3K56ac antibody (pAb)	ChIP, ChIP-Seq, DB, WB	100 µl	39281	¥70,000	→ ¥56,000	¥52,500
Histone H3K4me1 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µl	39297	¥70,000	→ ¥56,000	¥52,500
Histone H3K4ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µl	39381	¥70,000	→ ¥56,000	¥52,500
Histone H3, C-terminal antibody (pAb)	ELISA, WB	100 µl	39451	¥70,000	→ ¥56,000	¥52,500
Histone H2BK120ub1 antibody (mAb)	ChIP, WB	100 µg	39623	¥70,000	→ ¥56,000	¥52,500
Histone H4K20me3 antibody (mAb)	ChIP, DB, ICC, WB	100 µg	39671	¥70,000	→ ¥56,000	¥52,500
Histone H4R3me2a (asymmetric) antibody (pAb)	ChIP, DB, IF, WB	100 µl	39705	¥70,000	→ ¥56,000	¥52,500
Histone H2Av antibody (pAb)	ChIP, ChIP-Seq, WB	100 µl	39715	¥70,000	→ ¥56,000	¥52,500
Histone H4K20me1 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µg	39727	¥70,000	→ ¥56,000	¥52,500
Histone H3 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	100 µg	39763	¥70,000	→ ¥56,000	¥52,500
Histone H3K9me1 antibody (pAb)	ChIP, ChIP-Seq, DB, WB	100 µg	39887	¥70,000	→ ¥56,000	¥52,500
Histone H4K16ac antibody (pAb)	ChIP, DB, IHC, WB	100 µg	39929	¥70,000	→ ¥56,000	¥52,500
Histone H4ac (pan-acetyl) antibody (mAb)	ICC, IF, WB	100 µg	39967	¥70,000	→ ¥56,000	¥52,500
Histone H3K27me1 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µg	61015	¥70,000	→ ¥56,000	¥52,500
Histone H3K36me3 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	100 µg	61101	¥70,000	→ ¥56,000	¥52,500
Histone H3K9me0 antibody (mAb)	WB	100 µg	61399	¥70,000	→ ¥56,000	¥52,500
Histone H4 antibody (mAb)	ChIP, WB	100 µg	61521	¥70,000	→ ¥56,000	¥52,500
Histone H4K20ac antibody (mAb)	ChIP, DB	100 µg	61531	¥70,000	→ ¥56,000	¥52,500
Histone H4K20me2 antibody (mAb)	DB, WB	100 µg	61533	¥70,000	→ ¥56,000	¥52,500
クロマチン修飾 (200種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
BRD4 antibody (pAb)	ChIP, ChIP-Seq, IHC, WB	100 µl	39909	¥70,000	→ ¥56,000	¥52,500
CTCF antibody (pAb)	ChIP, ChIP-Seq, EMSA, ICC, IF, IHC, WB	100 µl	61311	¥70,000	→ ¥56,000	¥52,500
HDAC1 antibody (pAb)	ChIP, ChIP-Seq, WB	100 µg	40967	¥70,000	→ ¥56,000	¥52,500
HDAC2 antibody (mAb)	ChIP, ICC, IF, WB	100 µl	39533	¥70,000	→ ¥56,000	¥52,500
HDAC3 antibody (pAb)	ChIP, WB	100 µg	40968	¥70,000	→ ¥56,000	¥52,500
HDAC4 antibody (pAb)	ChIP, ChIP-Seq, WB	100 µg	40969	¥70,000	→ ¥56,000	¥52,500
HDAC5 antibody (pAb)	ChIP, IP, WB	100 µg	40970	¥70,000	→ ¥56,000	¥52,500
HDAC6 antibody (pAb)	ChIP, WB	100 µg	40971	¥70,000	→ ¥56,000	¥52,500
HDAC11 antibody (pAb)	WB	200 µl	39208	¥70,000	→ ¥56,000	¥52,500
HIRA antibody (mAb)	ICC, IF, WB	100 µg	39557	¥70,000	→ ¥56,000	¥52,500
HP1α antibody (pAb)	WB	200 µl	39295	¥70,001	→ ¥56,000	¥52,500
L3MBTL2 antibody (pAb)	IP, WB	100 µl	39569	¥70,000	→ ¥56,000	¥52,500
LSD1 / KDM1A antibody (pAb)	ChIP, WB	100 µl	61607	¥70,000	→ ¥56,000	¥52,500
MLL / HRX antibody (mAb)	ChIP, IP, WB	100 µg	39829	¥70,000	→ ¥56,000	¥52,500
MMSET / WHSC1 antibody (mAb)	ICC, IF, WB	100 µg	39879	¥70,000	→ ¥56,000	¥52,500
p300 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	100 µg	61401	¥70,000	→ ¥56,000	¥52,500
SMARCA2 / BRM antibody (mAb)	ChIP, ICC, IF, WB	100 µg	39805	¥70,000	→ ¥56,000	¥52,500

→裏面へ続く



リコンビナント抗体
AbFlex®も対象だよ!

一次抗体キャンペーン2本以下購入20% off 3本以上購入(組み合わせ自由)25% off

キャンペーン期間:2020年1月27日(月)~2020年3月31日(火) 受注分まで

DNAメチル化 (50種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
5-Carboxylcytosine (5-caC) antibody (pAb)	DB, ICC, IF, IHC, WB	100 µl	61225	¥70,000 →	¥56,000	¥52,500
5-Formylcytosine (5-fC) antibody (pAb)	DB, ICC, IF, IHC, WB	100 µl	61223	¥70,000 →	¥56,000	¥52,500
5-Hydroxymethylcytosine (5-hmC) antibody (pAb)	DB, FC, ICC, IF, IHC, MeDIP	100 µl	39769	¥70,000 →	¥56,000	¥52,500
5-Methylcytosine (5-mC) antibody (mAb)	DB, ELISA, ICC, IF, IHC, MeDIP, WB	50 µg	39649	¥70,000 →	¥56,000	¥52,500
BMI-1 antibody (mAb)	ChIP, ChIP-Seq, IF, IHC, WB	100 µg	39993	¥70,000 →	¥56,000	¥52,500
DNMT1 antibody (mAb)	ChIP, IHC, IP, WB	100 µg	39204	¥70,000 →	¥56,000	¥52,500
DNMT3A antibody (mAb)	ChIP, IHC, WB	100 µg	39206	¥70,000 →	¥56,000	¥52,500
DNMT3L antibody (pAb)	WB	100 µl	39907	¥70,000 →	¥56,000	¥52,500
EED antibody (mAb)	ChIP, ChIP-Seq, IHC	100 µg	61203	¥70,000 →	¥56,000	¥52,500
EZH2 antibody (pAb)	ChIP, ChIP-Seq, IHC, IP	100 µl	39901	¥70,000 →	¥56,000	¥52,500
Phc1 antibody (mAb)	ICC, IF, IP, WB	100 µg	39723	¥70,000 →	¥56,000	¥52,500
Phc2 antibody (mAb)	ChIP, ICC, IF, IP, WB	100 µg	39661	¥70,000 →	¥56,000	¥52,500
Ring1B antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	100 µg	39663	¥70,000 →	¥56,000	¥52,500
Suz12 antibody (pAb)	ChIP, ChIP-Seq, IF, WB	100 µl	39357	¥70,000 →	¥56,000	¥52,500
Tet1 antibody (pAb)	ChIP, ChIP-Seq, IHC	100 µl	61443	¥70,000 →	¥56,000	¥52,500
Tet2 antibody (mAb)	IP, WB	100 µg	61389	¥70,000 →	¥56,000	¥52,500
Tet3 antibody (pAb)	ChIP, WB	100 µl	61395	¥70,000 →	¥56,000	¥52,500
転写制御 (430種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
ATAD2 antibody (pAb)	IP, WB	100 µl	61369	¥70,000 →	¥56,000	¥52,500
CtIP antibody (mAb)	ICC, IF, IP, WB	100 µg	61141	¥70,000 →	¥56,000	¥52,500
HIF-1 alpha antibody (mAb)	IHC, WB	100 µg	61275	¥70,000 →	¥56,000	¥52,500
NF-1B2 antibody (pAb)	ChIP, EMSA, IF, WB	100 µg	39091	¥70,000 →	¥56,000	¥52,500
NF κB p65 antibody (pAb)	ChIP, ChIP-Seq, EMSA, WB	100 µl	39369	¥70,000 →	¥56,000	¥52,500
PLZF antibody (mAb)	ChIP, ICC, IF, IP, WB	100 µg	39987	¥70,000 →	¥56,000	¥52,500
RNA pol II antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	200 µl	39097	¥70,000 →	¥56,000	¥52,500
RNA pol II CTD phospho Ser2 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	100 µg	61083	¥70,000 →	¥56,000	¥52,500
RNA pol II CTD phospho Ser5 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	100 µg	61085	¥70,000 →	¥56,000	¥52,500
RNA pol II CTD phospho Ser7 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	100 µg	61087	¥70,000 →	¥56,000	¥52,500
RNA pol II CTD phospho Tyr1 antibody (mAb)	ChIP, IP, WB	100 µg	61383	¥70,000 →	¥56,000	¥52,500
YY1 antibody (pAb)	ChIP-Seq, IHC, IP, WB	100 µl	61779	¥70,000 →	¥56,000	¥52,500
核内レセプター (20種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
Androgen Receptor antibody (pAb)	ChIP, ChIP-Seq, WB	100 µl	39781	¥70,000 →	¥56,000	¥52,500
RAR-γ antibody (mAb)	EMSA, IHC, WB	100 µg	39973	¥70,000 →	¥56,000	¥52,500
DAX-1 / NR0B1 antibody (mAb)	ICC, IF, IHC, WB	100 µg	39983	¥70,000 →	¥56,000	¥52,500
RAR-α antibody (mAb)	EMSA, IHC, WB	100 µg	39971	¥70,000 →	¥56,000	¥52,500
RAR-β antibody (mAb)	EMSA, IHC, WB	100 µg	61027	¥70,000 →	¥56,000	¥52,500
Estrogen Receptor α antibody (mAb)	ChIP, IP, WB	100 µg	61035	¥70,000 →	¥56,000	¥52,500
LXR-α antibody (pAb)	ChIP, ChIP-Seq, WB	100 µl	61175	¥70,000 →	¥56,000	¥52,500
LXR-β antibody (pAb)	ChIP, ChIP-Seq, WB	100 µl	61177	¥70,000 →	¥56,000	¥52,500
RNAプロセッシング (10種以上)	適用	包装	商品コード	価格(税別)	キャンペーン価格(税別) 2本以下購入 20% off	3本以上購入 25% off
Ago1 antibody (mAb)	IP, WB	100 µg	61071	¥70,000 →	¥56,000	¥52,500
Ago1/2/3 antibody (mAb)	ChIP, ICC, IF, IHC, IP, WB	100 µg	39937	¥70,000 →	¥56,000	¥52,500
Ago2 antibody (mAb)	ICC, IF, IP, WB	100 µg	39853	¥70,000 →	¥56,000	¥52,500
Ago3 antibody (mAb)	ChIP, ICC, IF, IP	100 µg	39787	¥70,000 →	¥56,000	¥52,500
Ago4 antibody (mAb)	IP, WB	100 µg	39855	¥70,000 →	¥56,000	¥52,500
N6-Methyladenosine (m6A) antibody (pAb)	DB	100 µg	61495	¥70,000 →	¥56,000	¥52,500

略語:ChIP: Chromatin IP, ChIP-Seq: ChIP Sequencing, DB: Dot Blot, ELISA: Enzyme-Linked Immunosorbent Assay, EMSA: Electrophoretic Mobility Shift Assay, FACS: Fluorescence-Activated Cell Sorting, ICC: Immunocytochemistry, IF: Immunofluorescence, IHC: Immunohistochemistry, IP: Immunoprecipitation, MeDIP: Methyl-DNA IP, WB: Western Blot, pAb: polyclonal antibody, mAb: monoclonal antibody

掲載されていない抗体も多数ございます。抗体全リストにつきましては弊社ウェブサイト

<https://www.activemotif.com/catalog/64/primary-antibodies>をご覧ください。

バルク包装をご希望の方はお気軽に japantech@activemotif.com までお問い合わせください。

検索は
こちら➡



アクティブ・モティブ株式会社

取扱店

www.activemotif.jp

〒162-0824 東京都新宿区相場町2-21

Tel: 03-5225-3638 Fax: 03-5261-8733

e-mail: japantech@activemotif.com