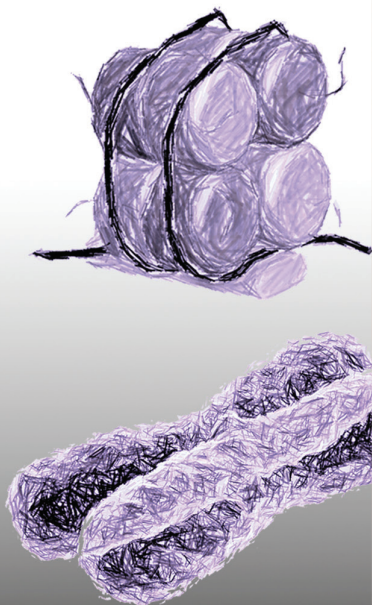


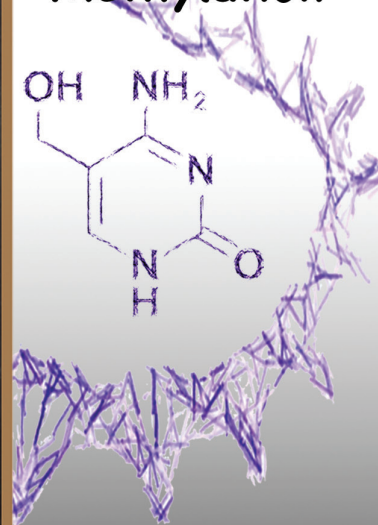
Histones



New to Epigenetics?
We can help!

ChIP Kits ChIP-SEQ
ChIP-validated Antibodies
DNA Methylation Variants
Histone Antibodies
Recombinant Histones
Epigenetic Services

DNA Methylation



Receive a FREE ANTIBODY* from Active Motif

Buy 2 antibodies and get an additional antibody of your choice for free.

Simply reference the code "AB2016" when you place the order.

* Free antibody must be the lowest priced antibody of the three ordered. Offer expires December 31, 2016 and limited to a single use per order and cannot be used with other promotions or discounts.

** Please click [HERE](#) for a list of eligible countries.

Active Motif offers a wide range of antibodies to epigenetic modifications, chromatin modifying proteins and transcription factors. They have been validated in applications such as ChIP-Seq, western blotting and immunofluorescence and many are available in a convenient sample size. For a complete list, please visit us at www.activemotif.com/abs.

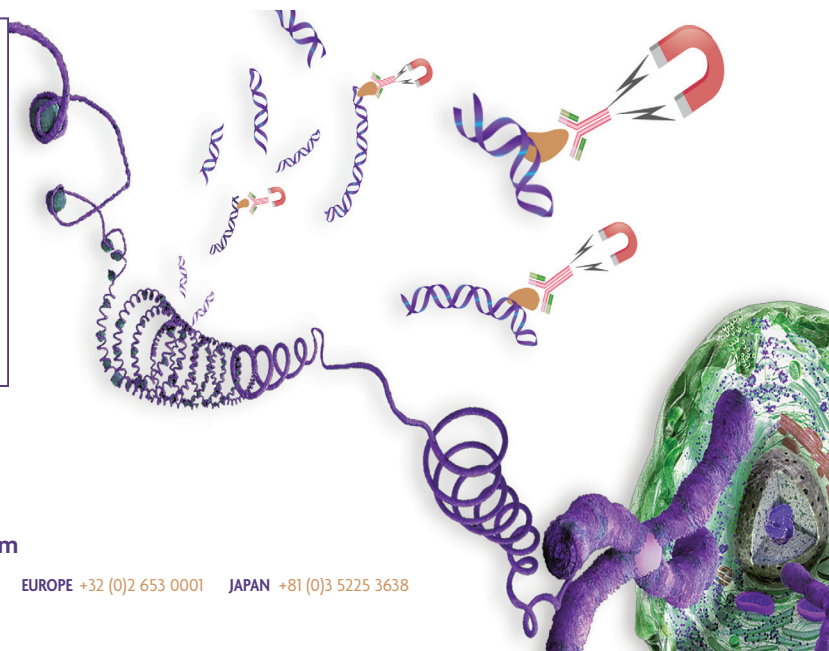
Simply reference the code "AB2016" when you place the order to orders@activemotif.com or contact your local **Key Account Manager** for more information. This offer is valid through December 31, 2016 and limited to a single use per order and cannot be used with other promotions or discounts.

Contact:

Active Motif

Email: orders@activemotif.com

877 222 9543



ACTIVE MOTIF®

Enabling Epigenetics Research

www.activemotif.com

NORTH AMERICA 877 222 9543 EUROPE +32 (0)2 653 0001 JAPAN +81 (0)3 5225 3638