

AMPIVIEW[®] Circ018 (AS) Dig RNA Probes (Mouse) Set

A digoxigenin-labeled RNA probe for the *in situ* hybridization (ISH) detection IncRNA

Circ018 in mouse or rat cells and tissues
AMPIVIEW[®] Circ018 (AS) Dig RNA Probes (mouse) are digoxigenin-labeled RNA probes targeting Circ018 circular RNA for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

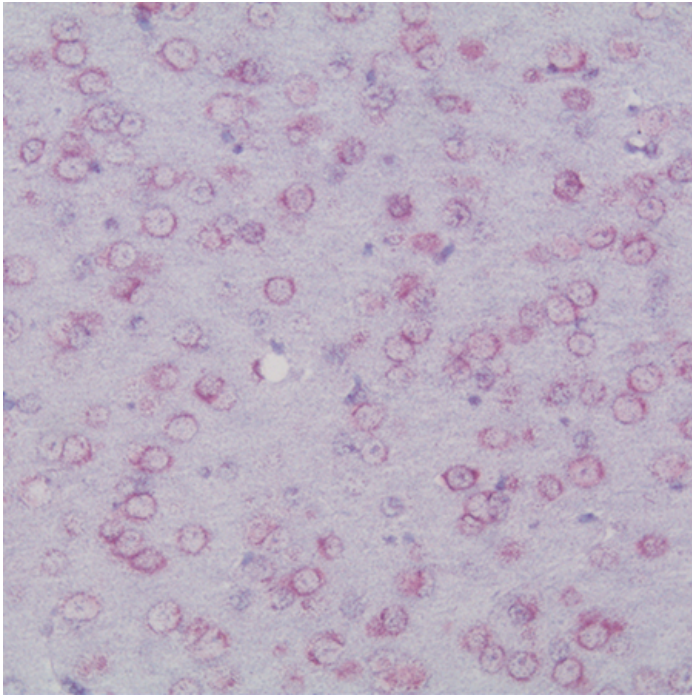
Ordering Information

[Order Online »](#)

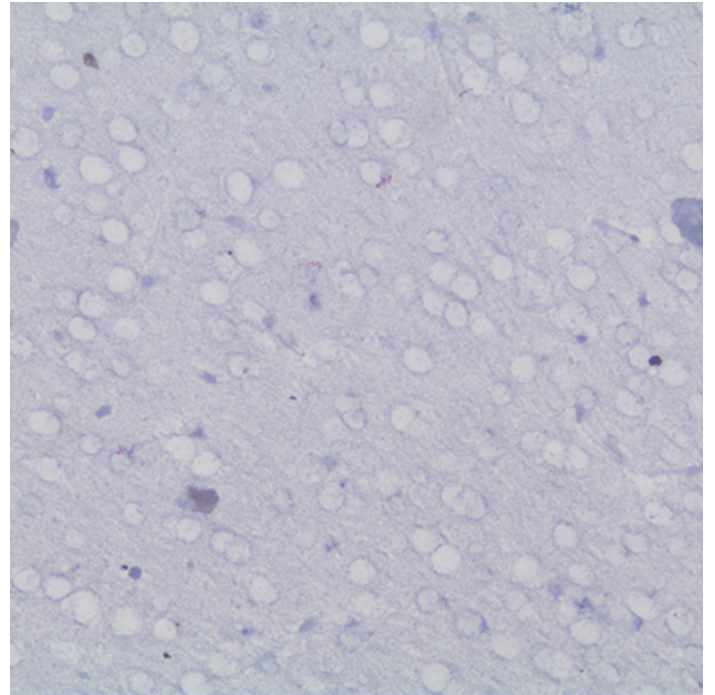
ENZ-GEN180-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Circ018 (red) detection in mouse brain tissue specimen with AMPIVIEW[®],_φ Circ018 (AS) Dig RNA probes amplified with DIGX[®] rabbit anti-digoxigenin linker (ENZ-ABS303), detected with POLYVIEW[®] PLUS AP (anti-rabbit) (ENZ-ACC110) combined with HIGHDEF[®] Red AP Substrate/Chromogen and hematoxylin.



Circ018 detection in mouse brain tissue specimen with AMPIVIEW[®],_φ NSP-Dig RNA probes (negative control), amplified with DIGX[®] rabbit anti-digoxigenin linker (ENZ-ABS303), detected with POLYVIEW[®] PLUS AP (anti-rabbit) (ENZ-ACC110) combined with HIGHDEF[®] Red AP Substrate/Chromogen and hematoxylin.

Handling & Storage

Use/Stability	Aliquot and store at -20°C or -80°C. Under these conditions, products are stable until its expiration dates.
Handling	Allow contents to warm up to room temperature prior to use.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application ISH (in situ hybridization)

Application Notes AMPIVIEW[®] Circ018 (AS) Dig RNA probes have been designed to target mouse or rat circular DNA 018 in tissues and cells. Circ018 contains a closed loop structure and lack free ends. Circ018 has recently been discovered to play a critical role in the diagnosis of certain cancers. AMPIVIEW[®] Circ018 (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX[®] anti-digoxigenin linker and nanopolymer-based detection systems, POLYVIEW[®] PLUS, combined with HIGHDEF[®] chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents 2 x 1 mL AMPIVIEW[®] Circ018 (AS) Dig RNA Probes (2µg/mL)
1 x 2 mL AMPIVIEW[®] Hybridization Buffer (1X)

Technical Info / Product Notes Circ018 is a circular RNA that has been recently discovered to play a critical role in certain cancers.

AMPIVIEW[®] Circ018 (AS) Dig RNA Probes are optimized for detection of ubiquitin RNA or RNA/DNA in FFPE tissues or cells with DIGX[®] anti-digoxigenin linker, POLYVIEW[®] PLUS detection reagents combined with HIGHDEF[®] chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW[®] Hybridization Buffer (1X), included in this set.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)