# AMPIVIEW® CD19 (AS) Dig RNA Probes Set

A digoxigenin-labeled RNA probe for the *in* situ hybridization (ISH) detection of CD19

probes targeting CD19 nucleic acid (DNA/RNA or RNA) for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

# **Ordering Information**

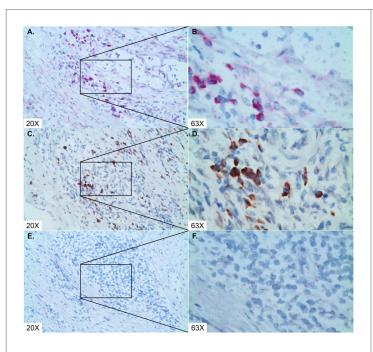
Order Online »

ENZ-GEN166-2000

2ml

Manuals, SDS & CofA

**View Online** »



CD19 was detected in colon, Burkitt's lymphoma with A.-D. AMPIVIEW® CD19 (AS) Dig RNA probes and E.-F. AMPIVIEW® NSP Dig RNA probes, amplified with DIGX® rabbit anti-digoxigenin linker and detected with A.-B. and E.-F. POLYVIEW® PLUS (anti-r) AP reagent combined with HIGHDEF® Red AP chromogen or C.D. POLYVIEW® PLUS (anti-r) HRP reagent combined with HIGHDEF® DAB chromogen/substrate and counterstained with HIGHDEF® hematoxylin.

## **Handling & Storage**

**Use/Stability** Upon receipt, aliquot and store at -20°C. For long-term storage, store at -80°C. These

products, under these conditions, are stable until the expiration date.

**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<sup>®</sup> CD19 (AS) Dig RNA probes have been

lymphocyte antigen CD19 in tissues and cells. CD19 is a transmembrane protein expressed in B cells and plasma cells. AMPIVIEW<sup>®</sup> CD19 (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to

be amplified with DIGX<sup>®</sup> anti-digoxigenin linker and detected with POLYVIEW<sup>®</sup> PLUS combined with

designed to detect the nucleic acid that encodes B-

HIGHDEF<sup>®</sup> chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents 2 x 1 mL AMPIVIEW® CD19 (AS) Dig RNA Probes

 $(2\mu g/mL)$ 

1 x 2 mL AMPIVIEW<sup>®</sup> Hybridization Buffer (1X)

**Technical Info / Product Notes**AMPIVIEW<sup>®</sup> CD19 (AS) Dig RNA Probes are optimized

for detection of CD19 RNA or RNA/DNA in FFPE tissues or cells with DIGX<sup>®</sup> anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> Hybridization

Buffer (1X), included in this set.



European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

m infoeu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com