

m1I antibody

Cat. No. C15410366

| | |
|--|--|
| Type: Polyclonal | Specificity: Human, mouse, other (wide range): positive |
| Size: 50 µg | Isotype: NA |
| Concentration: 2.1 µg/µl | Host: Rabbit |
| Lot No.: A2972-0010 | Purity: Protein A purified polyclonal antibody |
| Storage buffer: PBS containing 0.05% azide | Storage conditions: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles. |
| Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures. | |

Last Data Sheet Update: May 7, 2020

Description

Full name: **1-methylinosine polyclonal antibody**

Polyclonal antibody raised in rabbit against 1-methylinosine (m1I) conjugated to BSA.

Applications

| Applications | Suggested dilution | References |
|--------------|--------------------|------------|
| Dot Blot | 1:400 | Fig 1 |

Target Description

1-methylinosine (m1I) is a modified base which is mainly present in tRNAs.

Validation data

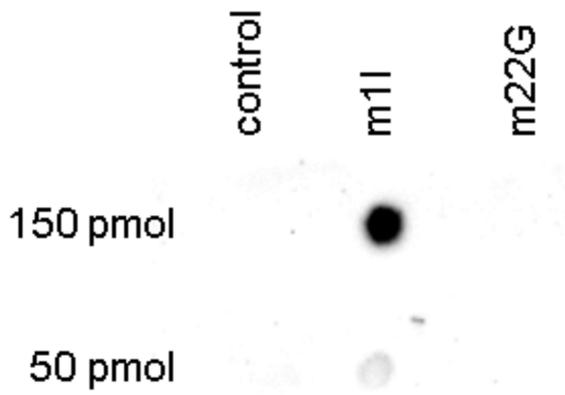


Figure 1. Dot blot analysis using the Diagenode antibody directed against m1I

To demonstrate the specificity of the Diagenode antibody against m1I (cat. No. C15410366), a Dot Blot analysis was performed using synthetic oligonucleotides containing different modified bases. 150 and 50 pmol of the respective oligo's were spotted on the membrane. The antibody was diluted of 1:400 in TBS-T containing 10 % skimmed milk and 1% BSA. Figure 1 shows a high specificity of the antibody for the oligonucleotide with the 1-methylinosine modification.