DATA SHEET





■ Parvovirus B19 VLP VP2

recombinant, Sf9 insect cells

Cat. No.	Amount
PR-1277	20 μg

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Purity: > 80 % (SDS-PAGE)

Form: liquid (Supplied in 16 mM sodium phosphate pH 7.4, 120 mM

NaCl and 20% glycerol)

Applications:

Western Blot with anti-Parvovirus B19 positive sera. Binds IgG-type human antibodies.

Description:

Recombinant Parvovirus B19 VLP VP2 produced in SF9 is a glycosylated, polypeptide chain having a calculated molecular mass of 60,839 Dalton.

The B19 virus, usually referred to as parvovirus B19 or sometimes as erythrovirus B19, belongs to the family of parvoviruses, genus erythrovirus. B19 virus causes a childhood rash named 'fifth disease' or 'erythema infectiosum' which is ordinarily known as 'slapped cheek syndrome'. Erythroviruses are members of the Parvoviridae family of small DNA viruses. The B19 virus is a non-enveloped, icosahedral virus which contains a single-stranded linear DNA genome. Parvovirus B19 is classified as erythrovirus because of its capability to invade red blood cell precursors in the bone marrow. This virus is mainly spread by infected respiratory droplets, though, blood-borne transmission has also been reported.

Selected References:

Beersma et al. (2005) Parvovirus B19 viral loads in relation to VP1 and VP2 antibody responses in diagnostic blood samples. J Clin Virol. **34**:71.

Manaresi et al. (2001) Differential IgM response to conformational and linear epitopes of parvovirus B19 VP1 and VP2 structural proteins. 64:67.