

**Chimeric Chagas Multiantigen His**

MACH  
recombinant, *E. coli*

Cat. No.	Amount
PR-1286	5 µg

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Purity:** > 90 % (SDS-PAGE)

**Form:** lyophilised (from 20mM Tris-HCl pH 8.5, 100mM sodium chloride, 20% trehalose and 0.1% sodium azide)

**Applications:**

Suitable for use in ELISA and Lateral flow.

**Description:**

Recombinant Chimeric Chagas Multiantigen (MACH) (*Trypanosoma cruzi*) produced in *E. coli*, is a polypeptide chain of 87 aa with epitopes PEP-2, TcD, TcE and SAPA. The protein is fused to a 6His-tag.

*Trypanosoma cruzi*, better known as *T. cruzi*, is a deadly parasite that causes Chagas' disease. This disease is a chronic infection, which primarily affects the heart and nervous system, causing severe neurological disorders, as well as swelling or denervation of nervous tissue in the heart, colon and esophagus. Chagas disease often goes undiagnosed due to close association of symptoms to heart disease and a variety of other disorders. The organism can circulate in the blood of infected patients for many years after infection, and can lead to transfusion-acquired infections in blood recipients from these infected donors. Contaminated blood transfusions are suspected to be the primary way in which the parasite has been transmitted to industrialized countries.

**Selected References:**

Umezawa *et al.* (2003) An improved serodiagnostic test for Chagas disease employing a mixture of *Trypanosoma cruzi* recombinant antigens. *Transfusion* 43:91.