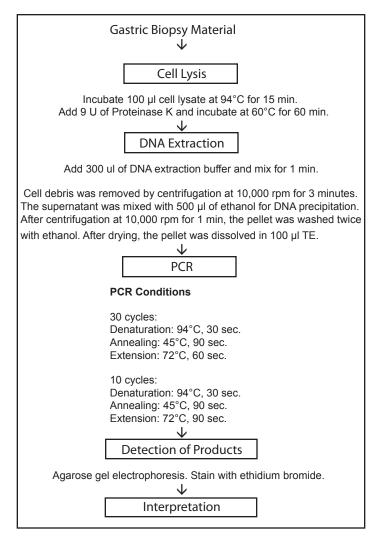
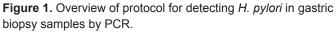
## Application: Comparison of the sensitivity of *TaKaRa Taq*<sup>™</sup> and *TaKaRa Ex Taq*<sup>®</sup> for detection of *Helicobacter pylori* in patient samples\*

*Helicobacter pylori* is the causative agent of gastritis and is associated with a variety of stomach disorders including gastric ulcers. Conventional detection of *H. pylori* in gastric biopsy specimens relies upon culture methods. PCR methods are now recognized as an effective alternative. PCR screening allows sensitive detection, and can provide results in much less time than culture-based methods.

The performance of *TaKaRa Taq*<sup>m</sup> and *TaKaRa Ex Taq*<sup> $\otimes$ </sup> in identifying *H. pylori* in patient samples was compared. Specifically, the polymerases were used to amplify the urease gene (HPUA) of *H. pylori* from gastric biopsy specimens from patients with gastritis. All patient samples were confirmed by culture methods to be positive for *H. pylori*-associated gastritis.





\*Data kindly provided by Dr. Kurokawa, Dr. Nukina, and Dr. Nakanishi (Public Health Research Institute of Kobe City)

# TakaRa

**TAKARA BIO INC**. 800-662-2566 Table 1. PCR Reaction Mixture Composition

Template DNA	10 µl
TaKaRa Ex Taq or Taq DNA Polymerase	2.5 U
10X Ex Taq Buffer or PCR Buffer	
dNTP Mixture	200 µM
Primers**	0.2 μM each

\*\*Primers to amplify HPUA

HPUA-1 (5'-GCCAATGGTAAATTAGTT 3': 304-321) HPUA-2 (5'-CTCCTTAATTGTTTTTAC 3': 714-697)

#### Results

Reactions performed with *TaKaRa Ex Taq* resulted in a distinct band for all of the biopsy samples, whereas the product was not amplified in all reactions when conventional *Taq* was used (Figure 2).

TaKaRa Ex Taq

Lane 5: H. pylori NCTC11637

56789

Lane 6: Gastric biopsy (1)

Lane 7: Gastric biopsy (2)

#### Taq

Lane 1: *H. pylori* NCTC11637 Lane 2: Gastric biopsy (1) Lane 3: Gastric biopsy (2) Lane 4: Gastric biopsy (3)

Lane 9: Marker

(3) Lane 8: Gastric biopsy (3)

**Figure 2.** Detection of *H. pylori* (HPUA) amplification products (410 bp). Three gastric biopsy samples (1, 2, 3) and *H. pylori* NCTC11637 (positive control, lane 1 and 5) were amplified with either *Taq* (lanes 1-4) or *Ex Taq* (lanes 5-8).

### Conclusion

The results indicate that *TaKaRa Ex Taq* was able to efficiently amplify the *H. pylori* HUPA target from patient specimens. These results suggest that *Ex Taq* minimizes the risk of falsenegative results that can occur using conventional *Taq* DNA polymerases.

Notice to Purchaser. Your use of these products and technologies is subject to compliance with any applicable licensing requirements described on the product's web page at http://www.clontech.com/takara. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements. The Takara logo is a trademark of TAKARA HOLDINGS, Kyoto, Japan. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.