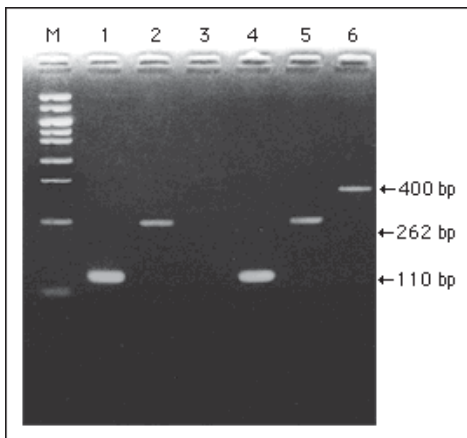


Application: Amplification of β -globin, K-ras 12, p53 (exon 5, exon 6) genes from 10-year-old paraffin sections

TaKaRa DEXPAT™ (DNA Extraction from Paraffin-embedded Tissue) (Cat.# 9091)

Genomic DNA was extracted from sections of 10-year-old paraffin-embedded colon cancer tissues using DEXPAT™ and used as a template for PCR to amplify fragments from the β -globin, K-ras 12 and p53 genes. To achieve accurate and highly sensitive PCR amplification, DNA extracted with DEXPAT™ is best used in conjunction with *TaKaRa Ex Taq™* (Cat. # RR001)

A. Amplification of β -globin gene fragments (110, 262 and 400 bp):

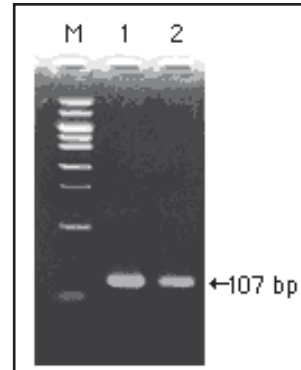


Lane M: pHY Marker (100 ng)
 Lane 1-3: Amplification products using 5 μ l of template DNA extracted using a conventional method
 Lane 4-5: Amplification products using 5 μ l of template DNA extracted using DEXPAT™
 PCR volume: 50 μ l
 Polymerase: *TaKaRa Ex Taq™*

Thermal Cycling Conditions:

94°C, 30 sec. }
 54°C, 60 sec. } 35 cycles
 72°C, 60 sec. }
 ↓
 72°C, 5 min.: 1 cycle

B. Amplification of K-ras gene (107 bp)

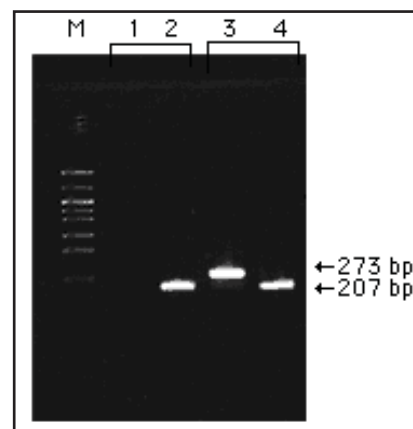


Lane M: pHY Marker (100 ng)
 Lane 1-3: Amplification products using 5 μ l of template DNA extracted using a conventional method
 PCR volume: 50 μ l
 Polymerase: *TaKaRa Ex Taq™*

Thermal Cycling Conditions:

94°C, 30 sec. }
 54°C, 60 sec. } 35 cycles
 72°C, 60 sec. }
 ↓
 72°C, 5 min.: 1 cycle

C. Amplification of p53 gene fragments from exons 5 (273 bp) and 6 (207 bp):



Lane M: pHY Marker (100 ng)
 Lane 1 & 2: Amplification products from DNA extracted using a conventional method
 Lane 3 & 4: Amplification products from DNA extracted using DEXPAT™
 PCR volume: 50 μ l
 Polymerase: *TaKaRa Ex Taq™*

continued ...

Application: Amplification of β -globin, K-ras 12, p53 (exon 5, exon 6) genes from 10-year-old paraffin sections

TaKaRa DEXPAT™ (DNA Extraction from Paraffin-embedded Tissue) (Cat.# 9091)

(continued)

1st PCR thermal cycling conditions:

Template: 5 μ l of template DNA extracted using DEXPAT™ or a conventional method
PCR: 50 μ l
Polymerase: TaKaRa Ex Taq™

94°C, 1 min..
60°C, 1 min.
72°C, 2 min. } 30 cycles
↓
72°C, 5 min.: 1 cycle

2nd PCR thermal cycling conditions:

Template: 1 μ l of 1st PCR product
PCR: 50 μ l
Polymerase: TaKaRa Ex Taq™

94°C, 30 sec.
60°C, 30 sec.
72°C, 60 sec. } 25 cycles
↓
72°C, 5 min.: 1 cycle