

PRODUCT NAME	
Arabidopsis FLC-Intron1 Primer Pair	
Other names: AGAMOUS-LIKE 25, AT5G10140, AGL25, FLC, FLF, FLOWERING LOCUS C, FLOWERING LOCUS F, REDUCED STEM BRANCHING 6, RSB6	
Cat. No: C17040001	Format: 50 µl
	Format: 500 µl

Description: This primer pair specifically amplifies a genomic region from intron 1 of the Arabidopsis FLC gene. These Primers are optimized for routine SYBR® Green Real-Time qPCR assay following ChIP

Amplicon length: 84 base pairs

Amplified region: chr5: 3,175,656-3,175,739 (UCSC Jan 2004)

Specificity: *Arabidopsis thaliana*

Format: 10 µM solution in MilliQ water (5 µM of each primer)

Storage: For long storage, store at -20°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: November 25, 2014

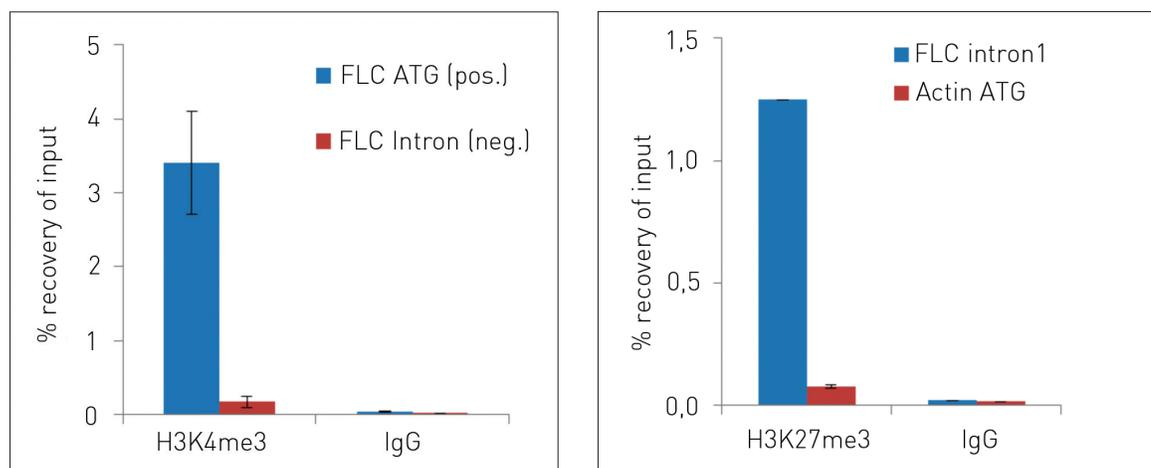


Figure 1. ChIP was performed on *Arabidopsis thaliana* (Col-0) seedlings 14 days after germination (dag) using our premium anti-H3K4me3 ChIP-seq grade antibody (#C15410003), anti-H3K27me3 ChIP-seq grade antibody (#C15410195) and rabbit IgG (#C15410206). Data shown is derived from biological replicates. Sheared chromatin from 0.2 g (fw) of seedlings was used per IP. Quantitative PCR was performed with the control FLC-ATG, FLC-Intron1 and Actin ATG primer sets from the kit. The recovery is expressed as % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).